

PRODUCTS AND SERVICES

Target list

Kinase Products

P.1-11 Kinase Products

Biochemical Assays

P.12 "QuickScout Screening Assist™ Kits" Kinase Protein Assay Kits

P.13 "QuickScout Custom Profiling & Panel Profiling Series" Targets

P.14 "QuickScout Custom Profiling Series" Preincubation Targets

Cell-Based Assays

P.15-16 NanoBRET™ TE Intracellular Kinase Cell-Based Assay Service Targets

P.17 Detection of Protein-Protein Interactions ~ProbeX™~ Stable Cell Lines

Kinase Products

Tyrosine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
ABL(ABL1)	HIS	ABL(ABL1)	08-001	Full-length	2-1130	NP_005148.2	N-terminal His	Insect (sf21)
	BTN	BTN-ABL(ABL1)	08-401-20N	Full-length	2-1130	NP_005148.2	N-terminal DYKDDDDK	Insect (sf21)
ABL(ABL1) [E255K]	HIS	ABL(ABL1)[E255K]	08-094	Full-length	2-1130	NP_005148.2	N-terminal His	Insect (sf21)
ABL(ABL1) [T315I]	HIS	ABL(ABL1)[T315I]	08-093	Full-length	2-1130	NP_005148.2	N-terminal His	Insect (sf21)
	BTN	BTN-ABL(ABL1)[T315I]	08-493-20N	Full-length	2-1130	NP_005148.2	N-terminal DYKDDDDK	Insect (sf21)
ACK(TNK2)	GST	ACK(TNK2)	08-196	Catalytic domain	110-476	NP_005772.3	N-terminal GST	Insect (sf21)
ALK	GST	ALK	08-518	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK	08-405-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [C1156Y]	GST	ALK[C1156Y]	08-530	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[C1156Y]	08-730-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [D1203N]	GST	ALK[D1203N]	08-583	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
ALK [F1174L]	GST	ALK[F1174L]	08-519	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[F1174L]	08-719-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [F1174L/G1202R]	GST	ALK[F1174L/G1202R]	08-568	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[F1174L/G1202R]	08-768-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [G1202R]	GST	ALK[G1202R]	08-544	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[G1202R]	08-744-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [G1269A]	GST	ALK[G1269A]	08-537	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[G1269A]	08-737-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [I1171N]	GST	ALK[I1171N]	08-566	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[I1171N]	08-766-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [I1171N/L1198F]	GST	ALK[I1171N/L1198F]	08-567	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[I1171N/L1198F]	08-767-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [I1171S]	GST	ALK[I1171S]	08-569	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[I1171S]	08-769-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [I1171T]	GST	ALK[I1171T]	08-590	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[I1171T]	08-790-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [L1196M]	GST	ALK[L1196M]	08-529	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[L1196M]	08-429-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [L1196M/G1202R]	GST	ALK[L1196M/G1202R]	08-556	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[L1196M/G1202R]	08-756-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [L1198F]	GST	ALK[L1198F]	08-571	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[L1198F]	08-771-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [L1198F/G1202R]	GST	ALK[L1198F/G1202R]	08-584	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
ALK [R1275Q]	GST	ALK[R1275Q]	08-520	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[R1275Q]	08-720-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ALK [T1151_L1152insT]	GST	ALK[T1151_L1152insT]	08-539	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
ALK [V1180L]	GST	ALK[V1180L]	08-570	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ALK[V1180L]	08-770-20N	Cytoplasmic domain	1058-1620	NP_004295.2	N-terminal DYKDDDDK	Insect (sf21)
ARG(ABL2)	GST	ARG(ABL2)	08-102	Full-length	2-1161	NP_001161708.1	N-terminal GST	Insect (sf21)
	GST	AXL	08-107	Cytoplasmic domain	464-885	NP_001690.2	N-terminal GST	Insect (sf21)
AXL	BTN	BTN-AXL	08-407-20N	Cytoplasmic domain	464-885	NP_001690.2	N-terminal DYKDDDDK	Insect (sf21)
BLK	GST	BLK	08-164	Full-length	1-505	NP_001706.2	N-terminal GST	Insect (sf21)
BMX	GST	BMX	08-179	Full-length	1-675	NP_001712.1	N-terminal GST	Insect (sf21)
	BTN	BTN-BMX	08-479-20N	Full-length	1-675	NP_001712.1	N-terminal DYKDDDDK	Insect (sf21)
BRK(PTK6)	GST	BRK(PTK6)	08-165	Full-length	2-451	NP_005966.1	N-terminal GST	Insect (sf21)
	BTN	BTN-BRK(PTK6)	08-465-20N	Full-length	2-451	NP_005966.1	N-terminal DYKDDDDK	Insect (sf21)
BTN-BRK(PTK6)[non-activated]	BTN	BTN-BRK(PTK6)[non-activated]	08-465-23N	Full-length	2-451	NP_005966.1	N-terminal DYKDDDDK	Insect (sf21)
BTK	GST	BTK	08-180	Full-length	2-659	NP_000052.1	N-terminal GST	Insect (sf21)
	BTN	BTN-BTK	08-480-20N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
BTK [A428D]	BTN	BTN-BTK[A428D]	08-480-23N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-BTK[A428D]	08-704-20N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
BTK [C481S]	GST	BTK[C481S]	08-547	Full-length	2-659	NP_000052.1	N-terminal GST	Insect (sf21)
	BTN	BTN-BTK[C481S]	08-417-20N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
BTK [M437R]	BTN	BTN-BTK[C481S][non-activated]	08-417-23N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	BTK[M437R]	08-805	Full-length	2-659	NP_000052.1	N-terminal GST	Insect (sf21)
BTK [M437R]	BTN	BTN-BTK[M437R]	08-705-20N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-BTK[M437R][non-activated]	08-705-23N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
BTK [T316A]	BTN	BTN-BTK[T316A]	08-418-20N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-BTK[T316A][non-activated]	08-418-23N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
BTK [T474I]	BTN	BTN-BTK[T474I]	08-419-20N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-BTK[T474I][non-activated]	08-419-23N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
BTK [T474S]	BTN	BTN-BTK[T474S]	08-420-20N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-BTK[T474S][non-activated]	08-420-23N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
BTK [V416L]	BTN	BTN-BTK[V416L]	08-706-20N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-BTK[V416L][non-activated]	08-706-23N	Full-length	2-659	NP_000052.1	N-terminal DYKDDDDK	Insect (sf21)
CSK	GST	CSK	08-111	Full-length	1-450	NP_004374.1	N-terminal GST	Insect (sf21)
DDR1	GST	DDR1	08-113	Cytoplasmic domain	444-876	NP_001945.3	N-terminal GST	Insect (sf21)
	BTN	BTN-DDR1	08-432-20N	Cytoplasmic domain	444-876	NP_001945.3	N-terminal DYKDDDDK	Insect (sf21)
DDR1(SRC treated)	GST	DDR1(SRC treated)	08-132	Cytoplasmic domain	444-876	NP_001945.3	N-terminal GST	Insect (sf21)
	BTN	BTN-DDR1(SRC treated)	08-732-20N	Cytoplasmic domain	444-876	NP_001945.3	N-terminal DYKDDDDK	Insect (sf21)
DDR2	GST	DDR2	08-114	Cytoplasmic domain	422-855	NP_006173.2	N-terminal GST	Insect (sf21)
	BTN	BTN-DDR2	08-414-20N	Cytoplasmic domain	422-855	NP_006173.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR	GST	EGFR	08-115	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR	08-415-20N	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [C797S]	GST	EGFR[C797S]	08-561	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[C797S]	08-761-20N	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [C797S/L858R]	GST	EGFR[C797S/L858R]	08-563	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[C797S/L858R]	08-763-20N	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [d746-750]	GST	EGFR[d746-750]	08-527	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[d746-750]	08-427-20N	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)

Kinase Products

Tyrosine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
EGFR [d746-750/C797S]	GST	EGFR[d746-750/C797S]	08-564	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[d746-750/C797S]	08-764-20N	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [d746-750/T790M]	GST	EGFR[d746-750/T790M]	08-528	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[d746-750/T790M]	08-428-20N	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [d746-750/T790M/C797S]	GST	EGFR[d746-750/T790M/C797S]	08-565	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [d746-750/T790M/C797S/L858R]	GST	EGFR[d746-750/T790M/C797S/L858R]	08-811	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[d746-750/T790M/C797S/L858R]	08-711-20N	Cytoplasmic domain	669-745, 751-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [D770_N771insNPG]	GST	EGFR[D770_N771insNPG]	08-553	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [D770_N771insNPG/T790M]	GST	EGFR[D770_N771insNPG/T790M]	08-558	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [L718M/T790M/C797S/L858R]	GST	EGFR[L718M/T790M/C797S/L858R]	08-801	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [L718Q/T790M/C797S/L858R]	GST	EGFR[L718Q/T790M/C797S/L858R]	08-802	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [L718V/T790M/C797S/L858R]	GST	EGFR[L718V/T790M/C797S/L858R]	08-803	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [L747P]	GST	EGFR[L747P]	08-588	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [L792H]	GST	EGFR[L792H]	08-562	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [L858R]	GST	EGFR[L858R]	08-502	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[L858R]	08-402-20N	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [L861Q]	GST	EGFR[L861Q]	08-513	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [N771_P772insH]	GST	EGFR[N771_P772insH]	08-149	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [N771GY]	GST	EGFR[N771GY]	08-110	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [S768I]	GST	EGFR[S768I]	08-589	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
EGFR [T790M]	GST	EGFR[T790M]	08-194	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[T790M]	08-494-20N	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [T790M/L858R]	GST	EGFR[T790M/L858R]	08-510	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[T790M/L858R]	08-410-20N	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EGFR [T790M/C797S/L858R]	GST	EGFR[T790M/C797S/L858R]	08-559	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EGFR[T790M/C797S/L858R]	08-759-20N	Cytoplasmic domain	669-1210	NP_005219.2	N-terminal DYKDDDDK	Insect (sf21)
EML4-ALK	GST	EML4-ALK	08-516	Fused	1-1059	BAF73611.1	N-terminal GST	Insect (sf21)
EPHA1	GST	EPHA1	08-119	Cytoplasmic domain	586-976	NP_005223.4	N-terminal GST	Insect (sf21)
EPHA2	GST	EPHA2	08-121	Cytoplasmic domain	572-976	NP_004422.2	N-terminal GST	Insect (sf21)
	BTN	BTN-EPHA2	08-421-20N	Cytoplasmic domain	572-976	NP_004422.2	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-EPHA2[non-activated]	08-421-23N	Cytoplasmic domain	572-976	NP_004422.2	N-terminal DYKDDDDK	Insect (sf21)
EPHA3	GST	EPHA3	08-122	Cytoplasmic domain	579-983	NP_005224.2	N-terminal GST	Insect (sf21)
EPHA4	GST	EPHA4	08-123	Cytoplasmic domain	586-986	NP_004429.1	N-terminal GST	Insect (sf21)
EPHA5	GST	EPHA5	08-124	Catalytic domain	662-948	NP_004430.4	N-terminal GST	Insect (sf21)
EPHA6	GST	EPHA6	08-125	Cytoplasmic domain	683-1130	NP_001073917.2	N-terminal GST	Insect (sf21)
EPHA7	GST	EPHA7	08-126	Cytoplasmic domain	595-998	NP_004431.1	N-terminal GST	Insect (sf21)
EPHA8	GST	EPHA8	08-127	Catalytic domain	571-924	NP_065387.1	N-terminal GST	Insect (sf21)
EPHB1	GST	EPHB1	08-128	Cytoplasmic domain	578-984	NP_004432.1	N-terminal GST	Insect (sf21)
EPHB2	GST	EPHB2	08-129	Cytoplasmic domain	581-987	NP_004433.2	N-terminal GST	Insect (sf21)
EPHB3	GST	EPHB3	08-130	Cytoplasmic domain	596-998	NP_004434.2	N-terminal GST	Insect (sf21)
EPHB4	GST	EPHB4	08-131	Cytoplasmic domain	577-987	NP_004435.3	N-terminal GST	Insect (sf21)
FAK(PTK2)	GST	FAK(PTK2)	08-137	Truncated	376-1052	NP_722560.1	N-terminal GST	Insect (sf21)
	BTN	BTN-FAK(PTK2)	08-437-20N	Truncated	376-1052	NP_722560.1	N-terminal DYKDDDDK	Insect (sf21)
FER	GST	FER	08-139	Full-length	1-822	NP_005237.2	N-terminal GST	Insect (sf21)
FES	GST	FES	08-140	Full-length	1-413, 416-822	NP_001996.1	N-terminal GST	Insect (sf21)
FGFR1	GST	FGFR1	08-133	Cytoplasmic domain	398-822	NP_075598.2	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR1	08-435-20N	Cytoplasmic domain	398-822	NP_075598.2	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR1[non-activated]	08-435-23N	Cytoplasmic domain	398-822	NP_075598.2	N-terminal DYKDDDDK	Insect (sf21)
FGFR1 [V561M]	GST	FGFR1[V561M]	08-536	Cytoplasmic domain	398-822	NP_075598.2	N-terminal GST	Insect (sf21)
FGFR2	GST	FGFR2	08-134	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2	08-434-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[non-activated]	08-434-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [C491A]	GST	FGFR2[C491A]	08-576	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[C491A]	08-903-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[C491A][non-activated]	08-903-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [C491A/V564I]	GST	FGFR2[C491A/V564I]	08-578	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[C491A/V564I]	08-905-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[C491A/V564I][non-activated]	08-905-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [C491A/V564L]	GST	FGFR2[C491A/V564L]	08-579	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[C491A/V564L]	08-906-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[C491A/V564L][non-activated]	08-906-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [E565A]	GST	FGFR2[E565A]	08-807	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[E565A]	08-907-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[E565A][non-activated]	08-907-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [K659M]	GST	FGFR2[K659M]	08-817	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[K659M]	08-927-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[K659M][non-activated]	08-927-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [L617V]	GST	FGFR2[L617V]	08-816	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[L617V]	08-926-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[L617V][non-activated]	08-926-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [M537I]	GST	FGFR2[M537I]	08-812	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[M537I]	08-922-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[M537I][non-activated]	08-922-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [M537I/M538L]	GST	FGFR2[M537I/M538L]	08-814	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[M537I/M538L]	08-924-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[M537I/M538L][non-activated]	08-924-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [M538L]	GST	FGFR2[M538L]	08-813	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[M538L]	08-923-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[M538L][non-activated]	08-923-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)

Kinase Products

Tyrosine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
FGFR2 [N549D]	GST	FGFR2[N549D]	08-808	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[N549D]	08-908-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[N549D][non-activated]	08-908-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [N549H]	GST	FGFR2[N549H]	08-532	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[N549H]	08-901-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[N549H][non-activated]	08-901-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [N549K]	GST	FGFR2[N549K]	08-809	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[N549K]	08-909-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[N549K][non-activated]	08-909-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [N549K/V564F]	GST	FGFR2[N549K/V564F]	08-815	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[N549K/V564F]	08-925-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[N549K/V564F][non-activated]	08-925-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [V564F]	GST	FGFR2[V564F]	08-810	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[V564F]	08-910-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[V564F][non-activated]	08-910-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [V564I]	GST	FGFR2[V564I]	08-546	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[V564I]	08-902-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[V564I][non-activated]	08-902-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR2 [V564L]	GST	FGFR2[V564L]	08-577	Cytoplasmic domain	399-821	NP_000132.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR2[V564L]	08-904-20N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FGFR2[V564L][non-activated]	08-904-23N	Cytoplasmic domain	399-821	NP_000132.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR3	GST	FGFR3	08-135	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR3	08-433-20N	Cytoplasmic domain	436-806	NP_000133.1	N-terminal DYKDDDDK	Insect (sf21)
FGFR3 [E587Q]	GST	FGFR3[E587Q]	08-821	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [G697C]	GST	FGFR3[G697C]	08-542	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [K650E]	GST	FGFR3[K650E]	08-501	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [K650M]	GST	FGFR3[K650M]	08-199	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [L608V]	GST	FGFR3[L608V]	08-822	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [N540K]	GST	FGFR3[N540K]	08-818	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [V553L]	GST	FGFR3[V553L]	08-819	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [V553M]	GST	FGFR3[V553M]	08-820	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [V555L]	GST	FGFR3[V555L]	08-548	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR3 [V555M]	GST	FGFR3[V555M]	08-543	Cytoplasmic domain	436-806	NP_000133.1	N-terminal GST	Insect (sf21)
FGFR4	GST	FGFR4	08-136	Cytoplasmic domain	460-802	NP_002002.3	N-terminal GST	Insect (sf21)
	BTN	BTN-FGFR4	08-436-20N	Cytoplasmic domain	460-802	NP_002002.3	N-terminal DYKDDDDK	Insect (sf21)
FGFR4 [N535K]	GST	FGFR4[N535K]	08-524	Cytoplasmic domain	460-802	NP_002002.3	N-terminal GST	Insect (sf21)
FGFR4 [V550E]	GST	FGFR4[V550E]	08-525	Cytoplasmic domain	460-802	NP_002002.3	N-terminal GST	Insect (sf21)
FGFR4 [V550L]	GST	FGFR4[V550L]	08-526	Cytoplasmic domain	460-802	NP_002002.3	N-terminal GST	Insect (sf21)
FGR	GST	FGR	08-166	Full-length	1-529	NP_005239.1	N-terminal GST	Insect (sf21)
FLT1	GST	FLT1	08-189	Cytoplasmic domain	781-1338	NP_002010.2	N-terminal GST	Insect (sf21)
	GST	FLT3	08-154	Cytoplasmic domain	564-993	NP_004110.2	N-terminal GST	Insect (sf21)
FLT3	BTN	BTN-FLT3	08-454-21N	Cytoplasmic domain	564-993	NP_004110.2	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FLT3[non-activated]	08-454-23N	Cytoplasmic domain	564-993	NP_004110.2	N-terminal DYKDDDDK	Insect (sf21)
FLT4	GST	FLT4	08-190	Cytoplasmic domain	798-1298	NP_002011.2	N-terminal GST	Insect (sf21)
	GST	FMS(CSF1R)	08-155	Cytoplasmic domain	538-972	NP_005202.2	N-terminal GST	Insect (sf21)
FMS(CSF1R)	BTN	BTN-FMS(CSF1R)	08-455-21N	Cytoplasmic domain	538-972	NP_005202.2	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-FMS(CSF1R)[non-activated]	08-455-20N	Cytoplasmic domain	538-972	NP_005202.2	N-terminal DYKDDDDK	Insect (sf21)
FRK	GST	FRK	08-167	Catalytic domain	223-505	NP_002022.1	N-terminal GST	Insect (sf21)
FYN [isoform a]	GST	FYN[isoforma]	08-168	Full-length	1-537	NP_002028.1	N-terminal GST	Insect (sf21)
	BTN	BTN-FYN[isoforma]	08-468-20N	Full-length	1-537	NP_002028.1	N-terminal DYKDDDDK	Insect (sf21)
FYN [isoform b]	GST	FYN[isoformb]	08-531	Full-length	1-534	NP_694592.1	N-terminal GST	Insect (sf21)
HCK	GST	HCK	08-169	Truncated	25-526	NP_002101.2	N-terminal GST	Insect (sf21)
HER2(ERBB2)	HIS	HER2(ERBB2)	08-016	Cytoplasmic domain	676-1255	NP_004439.1	N-terminal His	Insect (sf21)
HER4(ERBB4)	GST	HER4(ERBB4)	08-118	Cytoplasmic domain	676-1308	NP_005226.1	N-terminal GST	Insect (sf21)
IGF1R	GST	[IGF1R]	08-141	Cytoplasmic domain	959-1367	NP_000866.1	N-terminal GST	Insect (sf21)
	BTN	BTN-IGF1R	08-441-20N	Cytoplasmic domain	959-1367	NP_000866.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-IGF1R[non-activated]	08-441-23N	Cytoplasmic domain	959-1367	NP_000866.1	N-terminal DYKDDDDK	Insect (sf21)
INSR	GST	INSR	08-142	Catalytic domain	1005-1310	NP_000199.2	N-terminal GST	Insect (sf21)
	BTN	BTN-INSR	08-442-20N	Catalytic domain	1005-1310	NP_000199.2	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-INSR[non-activated]	08-442-23N	Catalytic domain	1005-1310	NP_000199.2	N-terminal DYKDDDDK	Insect (sf21)
IRR(INSRR)	GST	IRR(INSRR)	08-143	Cytoplasmic domain	953-1297	NP_050303.1	N-terminal GST	Insect (sf21)
ITK	GST	ITK	08-181	Full-length	2-620	NP_005537.3	N-terminal GST	Insect (sf21)
	BTN	BTN-ITK	08-481-20N	Full-length	2-620	NP_005537.3	N-terminal DYKDDDDK	Insect (sf21)
JAK1	GST	JAK1	08-144	Catalytic domain	850-1154	NP_002218.2	N-terminal GST	Insect (sf21)
	BTN	BTN-JAK1	08-444-20N	Catalytic domain	850-1154	NP_002218.2	N-terminal DYKDDDDK	Insect (sf21)
JAK2	HIS	JAK2	08-045	Catalytic domain	826-1132	NP_004963.1	N-terminal His	Insect (sf21)
	BTN	BTN-JAK2	08-445-20N	Catalytic domain	826-1132	NP_004963.1	N-terminal DYKDDDDK	Insect (sf21)
JAK2(JH1 JH2)	GST	JAK2(JH1JH2)	08-514	Catalytic/pseudo-kinase domain	532-1132	NP_004963.1	N-terminal GST	Insect (sf21)
JAK2(JH1 JH2) [V617F]	GST	JAK2(JH1JH2)[V617F]	08-515	Catalytic/pseudo-kinase domain	532-1132	NP_004963.1	N-terminal GST	Insect (sf21)
JAK3	HIS	JAK3	08-046	Catalytic domain	795-1124	NP_000206.2	N-terminal His	Insect (sf21)
KDR	GST	KDR	08-191	Cytoplasmic domain	790-1356	NP_002244.1	N-terminal GST	Insect (sf21)
	BTN	BTN-KDR	08-491-20N	Cytoplasmic domain	790-1356	NP_002244.1	N-terminal DYKDDDDK	Insect (sf21)
KIT	GST	KIT	08-156	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)
	BTN	BTN-KIT	08-456-21N	Cytoplasmic domain	544-976	NP_000213	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-KIT[non-activated]	08-456-23N	Cytoplasmic domain	544-976	NP_000213	N-terminal DYKDDDDK	Insect (sf21)
KIT [D816E]	GST	KIT[D816E]	08-541	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)
KIT [D816V]	GST	KIT[D816V]	08-505	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)
KIT [D816Y]	GST	KIT[D816Y]	08-534	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)
KIT [T670I]	GST	KIT[T670I]	08-195	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)
KIT [T670I/D816V]	GST	KIT[T670I/D816V]	08-555	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)
KIT [V560G]	GST	KIT[V560G]	08-504	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)

Kinase Products

Tyrosine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
KIT [V560G/D816V]	GST	KIT[V560G/D816V]	08-535	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)
KIT [V654A/D816V]	GST	KIT[V654A/D816V]	08-554	Cytoplasmic domain	544-976	NP_000213.1	N-terminal GST	Insect (sf21)
LCK	GST	LCK	08-170	Full-length	1-509	NP_005347.3	N-terminal GST	Insect (sf21)
	BTN	BTN-LCK	08-470-20N	Full-length	1-509	NP_005347.3	N-terminal DYKDDDDK	Insect (sf21)
LTK	GST	LTK	08-106	Catalytic domain	498-796	NP_002335.2	N-terminal GST	Insect (sf21)
LTK [I565N]	GST	LTK[I565N]	08-593	Catalytic domain	498-796	NP_002335.2	N-terminal GST	Insect (sf21)
LTK [L590M]	GST	LTK[L590M]	08-594	Catalytic domain	498-796	NP_002335.2	N-terminal GST	Insect (sf21)
LTK [L592F]	GST	LTK[L592F]	08-595	Catalytic domain	498-796	NP_002335.2	N-terminal GST	Insect (sf21)
LTK [L650F]	GST	LTK[L650F]	08-597	Catalytic domain	498-796	NP_002335.2	N-terminal GST	Insect (sf21)
LYNa	GST	LYNa	08-171	Full-length	1-512	NP_002341.1	N-terminal GST	Insect (sf21)
	BTN	BTN-LYNa	08-471-20N	Full-length	1-512	NP_002341.1	N-terminal DYKDDDDK	Insect (sf21)
LYNb	GST	LYNb	08-172	Full-length	1-491	NP_001104567.1	N-terminal GST	Insect (sf21)
	GST	MER(MERTK)	08-108	Cytoplasmic domain	528-999	NP_006334.2	N-terminal GST	Insect (sf21)
MER(MERTK)	BTN	BTN-MER(MERTK)	08-408-20N	Cytoplasmic domain	528-999	NP_006334.2	N-terminal DYKDDDDK	Insect (sf21)
	GST	MET	08-151	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal GST	Insect (sf21)
MET	BTN	BTN-MET	08-451-20N	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal DYKDDDDK	Insect (sf21)
	GST	MET[D1228H]	08-540	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal GST	Insect (sf21)
MET [D1228H]	BTN	BTN-MET[D1228H]	08-440-20N	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal DYKDDDDK	Insect (sf21)
	GST	MET[D1228N]	08-120	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal GST	Insect (sf21)
MET [D1228N]	BTN	BTN-MET[D1228N]	08-760-20N	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal DYKDDDDK	Insect (sf21)
	GST	MET[M1250T]	08-545	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal GST	Insect (sf21)
MET [M1250T]	BTN	BTN-MET[M1250T]	08-745-20N	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal DYKDDDDK	Insect (sf21)
	GST	MET[Y1230C]	08-709-20N	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal DYKDDDDK	Insect (sf21)
MET [Y1230C]	GST	MET[Y1230H]	08-178	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal GST	Insect (sf21)
	BTN	BTN-MET[Y1230H]	08-478-20N	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal DYKDDDDK	Insect (sf21)
MET [Y1230H]	GST	MET[Y1235D]	08-198	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal GST	Insect (sf21)
	BTN	BTN-MET[Y1235D]	08-498-20N	Cytoplasmic domain	956-1390	NP_000236.2	N-terminal DYKDDDDK	Insect (sf21)
MET [Y1235D]	GST	MUSK	08-153	Catalytic domain	527-869	NP_005583.1	N-terminal GST	Insect (sf21)
	BTN	BTN-MUSK	08-453-20N	Catalytic domain	527-869	NP_005583.1	N-terminal DYKDDDDK	Insect (sf21)
MUSK	BTN	BTN-MUSK[non-activated]	08-453-23N	Catalytic domain	527-869	NP_005583.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	NPM1-ALK	08-517	Fused	1-680	BAA08343.1	N-terminal GST	Insect (sf21)
NPM1-ALK	GST	PDGFRα(PDGFRα)	08-157	Cytoplasmic domain	550-1089	NP_006197.1	N-terminal GST	Insect (sf21)
	BTN	BTN-PDGFRα(PDGFRα)[non-activated]	08-457-23N	Cytoplasmic domain	550-1089	NP_006197.1	N-terminal DYKDDDDK	Insect (sf21)
PDGFRα(PDGFRα) [D842V]	GST	PDGFRα(PDGFRα)[D842V]	08-506	Cytoplasmic domain	550-1089	NP_006197.1	N-terminal GST	Insect (sf21)
PDGFRα(PDGFRα) [T674I]	GST	PDGFRα(PDGFRα)[T674I]	08-503	Cytoplasmic domain	550-1089	NP_006197.1	N-terminal GST	Insect (sf21)
PDGFRα(PDGFRα) [V561D]	GST	PDGFRα(PDGFRα)[V561D]	08-507	Cytoplasmic domain	550-1089	NP_006197.1	N-terminal GST	Insect (sf21)
PDGFRβ(PDGFRβ)	GST	PDGFRβ(PDGFRβ)	08-158	Cytoplasmic domain	557-1106	NP_002600.1	N-terminal GST	Insect (sf21)
	BTN	BTN-PDGFRβ(PDGFRβ)	08-458-20N	Cytoplasmic domain	557-1106	NP_002600.1	N-terminal DYKDDDDK	Insect (sf21)
PDGFRβ(PDGFRβ)	BTN	BTN-PDGFRβ(PDGFRβ)[non-activated]	08-458-23N	Cytoplasmic domain	557-1106	NP_002600.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	PYK2(PTK2B)	08-138	Full-length	1-967	NP_775267.1	N-terminal GST	Insect (sf21)
PYK2(PTK2B)	BTN	BTN-PYK2(PTK2B)	08-438-20N	Full-length	1-967	NP_775267.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	RET	08-159	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
RET	BTN	BTN-RET	08-459-20N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-RET[non-activated]	08-459-23N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
RET [G691S]	GST	RET[G691S]	08-522	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
	BTN	BTN-RET[G691S]	08-722-20N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
RET [G810C]	BTN	BTN-RET[G691S][non-activated]	08-722-23N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	RET[G810C]	08-572	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
RET [G810C/M918T]	GST	RET[G810C/M918T]	08-580	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
RET [G810R]	GST	RET[G810R]	08-573	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
RET [G810S]	GST	RET[G810S]	08-574	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
RET [M918T]	GST	RET[M918T]	08-508	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
	BTN	BTN-RET[M918T]	08-708-20N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
RET [M918T]	BTN	BTN-RET[M918T][non-activated]	08-708-23N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	RET[S891A]	08-523	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
RET [S891A]	BTN	BTN-RET[S891A]	08-723-20N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-RET[S891A][non-activated]	08-723-23N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
RET [V804M]	GST	RET[V804M]	08-560	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
	BTN	BTN-RET[V804M]	08-460-20N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
RET [V804M]	BTN	BTN-RET[V804M][non-activated]	08-460-23N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	RET[Y791F]	08-521	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal GST	Insect (sf21)
RET [Y791F]	BTN	BTN-RET[Y791F]	08-721-20N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-RET[Y791F][non-activated]	08-721-23N	Cytoplasmic domain	658-1114	NP_066124.1	N-terminal DYKDDDDK	Insect (sf21)
RON(MST1R)	GST	RON(MST1R)	08-152	Cytoplasmic domain	979-1400	NP_002438.2	N-terminal GST	Insect (sf21)
ROS(ROS1)	GST	ROS(ROS1)	08-163	Cytoplasmic domain	1883-2347	NP_002935.2	N-terminal GST	Insect (sf21)
ROS(ROS1) [G2032R]	GST	ROS(ROS1)[G2032R]	08-585	Cytoplasmic domain	1883-2347	NP_002935.2	N-terminal GST	Insect (sf21)
ROS(ROS1) [G2032R/L2086F]	GST	ROS(ROS1)[G2032R/L2086F]	08-587	Cytoplasmic domain	1883-2347	NP_002935.2	N-terminal GST	Insect (sf21)
ROS(ROS1) [L2086F]	GST	ROS(ROS1)[L2086F]	08-586	Cytoplasmic domain	1883-2347	NP_002935.2	N-terminal GST	Insect (sf21)
SRC	GST	SRC	08-173	Full-length	1-536	NP_005408.1	N-terminal GST	Insect (sf21)
	BTN	BTN-SRC	08-473-20N	Full-length	1-536	NP_005408.1	N-terminal DYKDDDDK	Insect (sf21)
SRM(SRMS)	GST	SRM(SRMS)	08-174	Catalytic domain	215-488	NP_543013.1	N-terminal GST	Insect (sf21)
SYK	GST	SYK	08-176	Full-length	1-635	NP_003168.2	N-terminal GST	Insect (sf21)
	BTN	BTN-SYK	08-476-20N	Full-length	1-635	NP_003168.2	N-terminal DYKDDDDK	Insect (sf21)
TEC	BTN	BTN-SYK[non-activated]	08-476-23N	Full-length	1-635	NP_003168.2	N-terminal DYKDDDDK	Insect (sf21)
	GST	TEC	08-182	Catalytic domain	359-631	NP_003206.2	N-terminal GST	Insect (sf21)
TIE2(TEK)	BTN	BTN-TEC	08-482-20N	Catalytic domain	359-631	NP_003206.2	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-TEC[non-activated]	08-482-23N	Catalytic domain	359-631	NP_003206.2	N-terminal DYKDDDDK	Insect (sf21)
TIE2(TEK)	GST	TIE2(TEK)	08-185	Cytoplasmic domain	771-1124	NP_000450.3	N-terminal GST	Insect (sf21)
	BTN	BTN-TIE2(TEK)	08-485-20N	Cytoplasmic domain	771-1124	NP_000450.3	N-terminal DYKDDDDK	Insect (sf21)
TNK1	GST	TNK1	08-104	Catalytic domain	106-390	NP_003976.2	N-terminal GST	Insect (sf21)

Kinase Products

Serine / Threonine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
TRKA(NTRK1)	GST	TRKA(NTRK1)	08-186	Cytoplasmic domain	436-790	NP_001012331.1	N-terminal GST	Insect (sf21)
	BTN	BTN-TRKA(NTRK1)	08-486-20N	Cytoplasmic domain	436-790	NP_001012331.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-TRKA(NTRK1)[non-activated]	08-486-23N	Cytoplasmic domain	436-790	NP_001012331.1	N-terminal DYKDDDDK	Insect (sf21)
TRKB(NTRK2)	GST	TRKB(NTRK2)	08-187	Cytoplasmic domain	456-822	NP_001018074.1	N-terminal GST	Insect (sf21)
	BTN	BTN-TRKB(NTRK2)	08-487-20N	Cytoplasmic domain	456-822	NP_001018074.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-TRKB(NTRK2)[non-activated]	08-487-23N	Cytoplasmic domain	456-822	NP_001018074.1	N-terminal DYKDDDDK	Insect (sf21)
TRKC(NTRK3)	GST	TRKC(NTRK3)	08-197	Cytoplasmic domain	456-825	NP_002521.2	N-terminal GST	Insect (sf21)
	BTN	BTN-TRKC(NTRK3)	08-497-20N	Cytoplasmic domain	456-825	NP_002521.2	N-terminal DYKDDDDK	Insect (sf21)
TXK	GST	TXK	08-183	Catalytic domain	260-527	NP_003319.2	N-terminal GST	Insect (sf21)
	BTN	BTN-TXK	08-483-20N	Catalytic domain	260-527	NP_003319.2	N-terminal DYKDDDDK	Insect (sf21)
TYK2	GST	TYK2	08-147	Catalytic domain	871-1187	NP_003322.3	N-terminal GST	Insect (sf21)
	BTN	BTN-TYK2	08-447-20N	Catalytic domain	871-1187	NP_003322.3	N-terminal DYKDDDDK	Insect (sf21)
TYK2(JH1 JH2)	GST	TYK2(JH1 JH2)	08-148	Catalytic/pseudo-kinase domain	556-1187	NP_003322.3	N-terminal GST	Insect (sf21)
	BTN	BTN-TYK2(JH1 JH2)	08-448-20N	Catalytic/pseudo-kinase domain	556-1187	NP_003322.3	N-terminal DYKDDDDK	Insect (sf21)
TYRO3	GST	TYRO3	08-109	Cytoplasmic domain	453-890	NP_006284.2	N-terminal GST	Insect (sf21)
	BTN	BTN-TYRO3	08-409-20N	Cytoplasmic domain	453-890	NP_006284.2	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-TYRO3[non-activated]	08-409-23N	Cytoplasmic domain	453-890	NP_006284.2	N-terminal DYKDDDDK	Insect (sf21)
YES(YES1)	GST	YES(YES1)	08-175	Full-length	1-543	NP_005424.1	N-terminal GST	Insect (sf21)
	BTN	BTN-YES(YES1)	08-475-20N	Full-length	1-543	NP_005424.1	N-terminal DYKDDDDK	Insect (sf21)
YES(YES1) [T348I]	GST	YES(YES1)[T348I]	08-533	Full-length	1-543	NP_005424.1	N-terminal GST	Insect (sf21)
ZAP70	GST	ZAP70	08-177	Full-length	1-619	NP_001070.2	N-terminal GST	Insect (sf21)

Serine / Threonine Kinases

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
ACTR2B(ACVR2B)	GST	ACTR2B(ACVR2B)	09-132	Cytoplasmic domain	161-512	NP_001097.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ACTR2B(ACVR2B)	09-432-20N	Cytoplasmic domain	161-512	NP_001097.2	N-terminal DYKDDDDK	Insect (sf21)
ACVR1B(ALK4)	GST	ACVR1B(ALK4)	09-135	Cytoplasmic domain	150-505	NP_004293.1	N-terminal GST	Insect (sf21)
	BTN	BTN-ACVR1B(ALK4)	09-435-20N	Cytoplasmic domain	150-505	NP_004293.1	N-terminal DYKDDDDK	Insect (sf21)
ACVR1C(ALK7)	GST	ACVR1C(ALK7)	09-150	Cytoplasmic domain	139-493	NP_060302.2	N-terminal GST	Insect (sf21)
ACVR2A	GST	ACVR2A	09-131	Cytoplasmic domain	162-513	NP_001607.1	N-terminal GST	Insect (sf21)
	BTN	BTN-ACVR2A	09-431-20N	Cytoplasmic domain	162-513	NP_001607.1	N-terminal DYKDDDDK	Insect (sf21)
ACVRL1(ALK1)	GST	ACVRL1(ALK1)	09-133	Truncated	139-503	NP_000011.2	N-terminal GST	Insect (sf21)
AKT1	GST	AKT1	01-101	Catalytic domain	104-480	NP_005154.2	N-terminal GST	Insect (sf21)
	BTN	BTN-AKT1	01-401-20N	Catalytic domain	104-480	NP_005154.2	N-terminal DYKDDDDK	Insect (sf21)
AKT2	GST	AKT2	01-102	Catalytic domain	120-481	NP_001617.1	N-terminal GST	Insect (sf21)
	BTN	BTN-AKT2	01-402-20N	Catalytic domain	120-481	NP_001617.1	N-terminal DYKDDDDK	Insect (sf21)
AKT3	GST	AKT3	01-103	Catalytic domain	108-479	NP_005456.1	N-terminal GST	Insect (sf21)
ALK2(ACVR1)	GST	ALK2(ACVR1)	09-134	Truncated	145-509	NP_001096.1	N-terminal GST	Insect (sf21)
ALK2(ACVR1) [Q207D]	GST	ALK2(ACVR1)[Q207D]	09-149	Truncated	145-509	NP_001096.1	N-terminal GST	Insect (sf21)
ALK2(ACVR1) [R206H]	GST	ALK2(ACVR1)[R206H]	09-148	Truncated	145-509	NP_001096.1	N-terminal GST	Insect (sf21)
AMPKα1/β1/γ1(PRKAA1/B1/G1)	GST	AMPKα1/β1/γ1(PRKAA1/B1/G1)	02-113	Full-length	1-550	NP_006242.5	N-terminal GST	Insect (sf21)
	BTN	BTN-AMPKα1/β1/γ1(PRKAA1/B1/G1)	02-413-20N	Full-length	1-550	NP_006242.5	N-terminal DYKDDDDK	Insect (sf21)
AMPKα1/β2/γ1(PRKAA1/B2/G1)	GST	AMPKα1/β2/γ1(PRKAA1/B2/G1)	02-147	Full-length	1-550	NP_006242.5	N-terminal GST	Insect (sf21)
	BTN	BTN-AMPKα1/β2/γ1(PRKAA1/B2/G1)	02-447-20N	Full-length	1-550	NP_006242.5	N-terminal DYKDDDDK	Insect (sf21)
AMPKα2/β1/γ1(PRKAA2/B1/G1)	GST	AMPKα2/β1/γ1(PRKAA2/B1/G1)	02-114	Full-length	1-552	NP_006243.2	N-terminal GST	Insect (sf21)
AMPKα2/β2/γ1(PRKAA2/B2/G1)	GST	AMPKα2/β2/γ1(PRKAA2/B2/G1)	02-148	Full-length	1-552	NP_006243.2	N-terminal GST	Insect (sf21)
AurA(AURKA)	GST	AurA(AURKA)	05-101	Full-length	1-403	NP_040835.1	N-terminal GST	Insect (sf21)
	BTN	BTN-AurA(AURKA)	05-401-20N	Full-length	1-403	NP_040835.1	N-terminal DYKDDDDK	Insect (sf21)
AurB(AURKB)/INCENP	GST	AurB(AURKB)/INCENP	05-102	Full-length	1-344	NP_004208.2	N-terminal GST	Insect (sf21)
	BTN	BTN-AurB(AURKB)/INCENP	05-402-21N	Full-length	1-344	NP_004208.2	N-terminal DYKDDDDK	Insect (sf21)
AurC(AURKC)	GST	AurC(AURKC)	05-103	Full-length	1-275	NP_003151.2	N-terminal GST	Insect (sf21)
BARK1(ADRBK1)	GST	BARK1(ADRBK1)	01-111	Full-length	1-689	NP_001610.2	N-terminal GST	Insect (sf21)
	BTN	BTN-BARK1(ADRBK1)	01-411-20N	Full-length	1-689	NP_001610.2	N-terminal DYKDDDDK	Insect (sf21)
BARK2(ADRBK2)	GST	BARK2(ADRBK2)	01-112	Full-length	1-688	NP_005151.2	N-terminal GST	Insect (sf21)
BMPRI1A	GST	BMPRI1A	09-137	Catalytic domain	187-532	NP_004320.2	N-terminal GST	Insect (sf21)
BMPRI1B	GST	BMPRI1B	09-138	Cytoplasmic domain	149-502	NP_001194.1	N-terminal GST	Insect (sf21)
BMPR2	GST	BMPR2	09-139	Cytoplasmic domain	174-1038	NP_001195.2	N-terminal GST	Insect (sf21)
BRAF	GST	BRAF	09-122	Catalytic domain	433-726	NP_004324.2	N-terminal GST	Insect (sf21)
	BTN	BTN-BRAF	09-422-20N	Catalytic domain	433-726	NP_004324.2	N-terminal DYKDDDDK	Insect (sf21)
BRAF [T599_V600insT]	GST	BRAF[T599_V600insT]	09-147	Catalytic domain	433-726	NP_004324.2	N-terminal GST	Insect (sf21)
BRAF [V600E]	GST	BRAF[V600E]	09-144	Catalytic domain	433-726	NP_004324.2	N-terminal GST	Insect (sf21)
BRSK1	GST	BRSK1	02-115	Full-length	1-778	NP_115806.1	N-terminal GST	Insect (sf21)
BRSK2	GST	BRSK2	02-116	Full-length	1-674	NP_001243558.1	N-terminal GST	Insect (sf21)
BUB1/BUB3	GST	BUB1/BUB3	05-187	Full-length	1-1085	NP_004327	N-terminal GST	Insect (sf21)
	BTN	BTN-BUB1/BUB3	05-487-20N	Full-length	1-1085	NP_004327	N-terminal DYKDDDDK	Insect (sf21)
BUBR1(BUB1B)	GST	BUBR1(BUB1B)	05-105	Full-length	1-1050	NP_001202.5	N-terminal GST	Insect (sf21)
CaMK1α(CAMK1)	GST	CaMK1α(CAMK1)	02-104	Full-length	1-370	NP_003647.1	N-terminal GST	Insect (sf21)
CaMK1β(PNCK)	GST	CaMK1β(PNCK)	02-105	Full-length	1-343	NP_001353906.1	N-terminal GST	Insect (sf21)
CaMK1δ(CAMK1D)	GST	CaMK1δ(CAMK1D)	02-106	Full-length	1-357	NP_065130.1	N-terminal GST	Insect (sf21)
CaMK2α(CAMK2A)	GST	CaMK2α(CAMK2A)	02-109	Full-length	1-478	NP_001196.1	N-terminal GST	Insect (sf21)
CaMK2β(CAMK2B)	GST	CaMK2β(CAMK2B)	02-110	Full-length	1-503	NP_001196.1(iso5)	N-terminal GST	Insect (sf21)
	BTN	BTN-CaMK2β(CAMK2B)	02-410-20N	Full-length	1-503	NP_001196.1(iso5)	N-terminal DYKDDDDK	Insect (sf21)
CaMK2γ(CAMK2G)	GST	CaMK2γ(CAMK2G)	02-112	Full-length	1-518	NP_001196.1	N-terminal GST	Insect (sf21)
CaMK2δ(CAMK2D)	GST	CaMK2δ(CAMK2D)	02-111	Full-length	1-478	NP_001196.1	N-terminal GST	Insect (sf21)
	BTN	BTN-CaMK2δ(CAMK2D)	02-411-20N	Full-length	1-478	NP_001196.1	N-terminal DYKDDDDK	Insect (sf21)
CaMK4	GST	CaMK4	02-108	Full-length	1-473	NP_001735.1	N-terminal GST	Insect (sf21)
CAMKK1	GST	CAMKK1	05-107	Full-length	1-505	NP_115670.1	N-terminal GST	Insect (sf21)
CAMKK2	GST	CAMKK2	05-108	Full-length	1-588	NP_006540.3(isoform-1)	N-terminal GST	Insect (sf21)
CDC7/ASK	GST	CDC7/ASK	05-109	Full-length	1-574	NP_003494.1	N-terminal GST	Insect (sf21)
CDK1(CDC2)	BTN	BTN-CDK1(CDC2)	04-480-20N	Full-length	1-297	NP_001777.1	N-terminal DYKDDDDK	Insect (sf21)

Kinase Products

Serine / Threonine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
CDK1(CDC2)/CycA2	GST	CDK1(CDC2)/CycA2	04-119	Full-length	1-297	NP_001777.1	N-terminal GST	Insect (sf21)
	BTN	BTN-CDK1(CDC2)/CycA2	04-419-20N	Full-length	1-297	NP_001777.1	N-terminal DYKDDDDK	Insect (sf21)
CDK1(CDC2)/CycB1	GST	CDK1(CDC2)/CycB1	04-102	Full-length	1-297	NP_001777.1	N-terminal GST	Insect (sf21)
	BTN	BTN-CDK1(CDC2)/CycB1	04-402-20N	Full-length	1-297	NP_001777.1	N-terminal DYKDDDDK	Insect (sf21)
CDK1(CDC2)/CycE1	GST	CDK1(CDC2)/CycE1	04-176	Full-length	1-297	NP_001777.1	N-terminal GST	Insect (sf21)
CDK2	BTN	BTN-CDK2	04-422-20N	Full-length	1-298	NP_001789.2	N-terminal DYKDDDDK	Insect (sf21)
CDK2/CycA2	GST	CDK2/CycA2	04-103	Full-length	1-298	NP_001789.2	N-terminal GST	Insect (sf21)
	BTN	BTN-CDK2/CycA2	04-403-20N	Full-length	1-298	NP_001789.2	N-terminal DYKDDDDK	Insect (sf21)
CDK2/CycE1	GST	CDK2/CycE1	04-165	Full-length	1-298	NP_001789.2	N-terminal GST	Insect (sf21)
	BTN	BTN-CDK2/CycE1	04-465-20N	Full-length	1-298	NP_001789.2	N-terminal DYKDDDDK	Insect (sf21)
CDK3	BTN	BTN-CDK3	04-481-20N	Full-length	1-305	NP_001249.1	N-terminal DYKDDDDK	Insect (sf21)
CDK3/CycC	GST	CDK3/CycC	04-174	Full-length	1-305	NP_001249.1	N-terminal GST	Insect (sf21)
	GST	CDK3/CycE1	04-104	Full-length	1-305	NP_001249.1	N-terminal GST	Insect (sf21)
CDK3/CycE1	BTN	BTN-CDK3/CycE1	04-404-20N	Full-length	1-305	NP_001249.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK4/CycD1	04-120	Full-length	1-303	NP_000066.1	N-terminal GST	Insect (sf21)
CDK4/CycD3	GST	CDK4/CycD3	04-105	Full-length	1-303	NP_000066.1	N-terminal GST	Insect (sf21)
	BTN	BTN-CDK4/CycD3	04-405-20N	Full-length	1-303	NP_000066.1	N-terminal DYKDDDDK	Insect (sf21)
CDK5	BTN	BTN-CDK5	04-493-20N	Full-length	1-292	NP_004926.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK5/p25	04-106	Full-length	1-292	NP_004926.1	N-terminal GST	Insect (sf21)
CDK5/p25	BTN	BTN-CDK5/p25	04-406-20N	Full-length	1-292	NP_004926.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK5/p35	04-192	Full-length	1-292	NP_004926.1	N-terminal GST	Insect (sf21)
CDK5/p35	BTN	BTN-CDK5/p35	04-492-20N	Full-length	1-292	NP_004926.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-CDK6	04-494-20N	Full-length	1-326	NP_001250.1	N-terminal DYKDDDDK	Insect (sf21)
CDK6	GST	CDK6/CycD1	04-114	Full-length	1-326	NP_001250.1	N-terminal GST	Insect (sf21)
CDK6/CycD3	GST	CDK6/CycD3	04-107	Full-length	1-326	NP_001250.1	N-terminal GST	Insect (sf21)
	BTN	BTN-CDK6/CycD3	04-407-20N	Full-length	1-326	NP_001250.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-CDK6/CycD3[non-activated]	04-407-23N	Full-length	1-326	NP_001250.1	N-terminal DYKDDDDK	Insect (sf21)
CDK7	BTN	BTN-CDK7	04-495-20N	Full-length	1-346	NP_001790.1	N-terminal DYKDDDDK	Insect (sf21)
CDK7/CycH/MAT1	GST	CDK7/CycH/MAT1	04-108	Full-length	1-346	NP_001790.1	N-terminal GST	Insect (sf21)
	BTN	BTN-CDK7/CycH/MAT1	04-408-20N	Full-length	1-346	NP_001790.1	N-terminal DYKDDDDK	Insect (sf21)
CDK8	BTN	BTN-CDK8	04-496-20N	Full-length	1-464	NP_001251.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK8/CycC	04-109	Full-length	1-464	NP_001251.1	N-terminal GST	Insect (sf21)
CDK8/CycC	BTN	BTN-CDK8/CycC	04-409-20N	Full-length	1-464	NP_001251.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK9/CycK	04-111	Full-length	1-372	NP_001252.1	N-terminal GST	Insect (sf21)
CDK9/CycK	BTN	BTN-CDK9/CycK	04-411-20N	Full-length	1-372	NP_001252.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK9/CycT1	04-110	Full-length	1-372	NP_001252.1	N-terminal GST	Insect (sf21)
CDK9/CycT1	BTN	BTN-CDK9/CycT1	04-410-20N	Full-length	1-372	NP_001252.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK9/CycT2	04-121	Full-length	1-372	NP_001252.1	N-terminal GST	Insect (sf21)
CDK9/CycT2	BTN	BTN-CDK9/CycT2	04-421-20N	Full-length	1-372	NP_001252.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK10(PISSLRE)/CycQ	04-171	Full-length	1-360	NP_443714.3	N-terminal GST	Insect (sf21)
CDK10(PISSLRE)/CycQ	BTN	BTN-CDK10(PISSLRE)/CycQ	04-471-20N	Full-length	1-360	NP_443714.3	N-terminal DYKDDDDK	Insect (sf21)
CDK12(CRKRS)(720-1490aa)/CycK	BTN	BTN-CDK12(CRKRS)(720-1490aa)/CycK	04-413-20N	Truncated	720-1490	NP_057591.2	N-terminal DYKDDDDK	Insect (sf21)
CDK13(CHED)(694-1512aa)/CycK	BTN	BTN-CDK13(CHED)(694-1512aa)/CycK	04-482-20N	Truncated	694-1512	NP_003709.3	N-terminal DYKDDDDK	Insect (sf21)
CDK13(CHED)/CycK	GST	CDK13(CHED)/CycK	04-123	Catalytic domain	694-1039	NP_003709.3	N-terminal GST, N-terminal His	Insect (sf21)
	BTN	BTN-CDK13(CHED)/CycK	04-483-20N	Catalytic domain	694-1039	NP_003709.3	N-terminal DYKDDDDK	Insect (sf21)
CDK19(CDC2L6)	BTN	BTN-CDK19(CDC2L6)	04-730-20N	Full-length	1-502	NP_055891.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CDK19(CDC2L6)/CycC	04-112	Full-length	1-502	NP_055891.1	N-terminal GST	Insect (sf21)
CDK19(CDC2L6)/CycC	BTN	BTN-CDK19(CDC2L6)/CycC	04-412-20N	Full-length	1-502	NP_055891.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CGK2(PRKG2)	01-143	Full-length	1-762	NP_006250.1	N-terminal GST	Insect (sf21)
CGK2(PRKG2)	GST	CHAK1(TRPM7)	10-111	Catalytic domain	1158-1865	NP_060142.3	N-terminal GST	Insect (sf21)
CHAK1(TRPM7)	GST	CHK1(CHEK1)	02-117	Full-length	1-476	NP_001265.2	N-terminal GST	Insect (sf21)
	BTN	BTN-CHK1(CHEK1)	02-417-20N	Full-length	1-476	NP_001265.2	N-terminal DYKDDDDK	Insect (sf21)
CHK1(CHEK1)	GST	CHK2(CHEK2)	02-162	Full-length	1-543	NP_009125.1	N-terminal GST	Insect (sf21)
CHK2(CHEK2)	GST	CK1α(CSNK1A1)	03-101	Full-length	1-337	NP_001883.4	N-terminal GST	Insect (sf21)
CK1α(CSNK1A1)	GST	CK1γ1(CSNK1G1)	03-105	Full-length	1-422	NP_071331.2	N-terminal GST	Insect (sf21)
CK1γ1(CSNK1G1)	GST	CK1γ2(CSNK1G2)	03-106	Full-length	1-415	NP_001310.3	N-terminal GST	Insect (sf21)
CK1γ2(CSNK1G2)	GST	CK1γ3(CSNK1G3)	03-107	Full-length	1-447	NP_004375.2	N-terminal GST	Insect (sf21)
CK1γ3(CSNK1G3)	GST	CK1δ(CSNK1D)	03-103	Catalytic domain	1-294	NP_001884.2	N-terminal GST	E. coli
CK1δ(CSNK1D)	GST	CK1ε(CSNK1E)	03-104	Catalytic domain	1-348	NP_001885.1	N-terminal GST	Insect (sf21)
CK1ε(CSNK1E)	GST	CK2α1/β(CSNK2A1/B)	05-184	Full-length(CK2α1)	1-391(CK2α1), 1-215(CK2β)	NP_001886.1(CK2α1), NP_001311.3(CK2β)	N-terminal GST(CK2α1), N-terminal His(CK2β)	Insect (sf21)
CK2α1/β(CSNK2A1/B)	GST/HIS	CK2α2/β(CSNK2A2/B)	05-185	Full-length(CK2α2)	1-350(CK2α2), 1-215(CK2β)	NP_001887.1(CK2α2), NP_001311.3(CK2β)	N-terminal His(CK2α2), N-terminal His(CK2β)	Insect (sf21)
CK2α2/β(CSNK2A2/B)	GST/HIS	CLK1	04-126	Catalytic domain	129-484	NP_004062.2	N-terminal GST	Insect (sf21)
CLK1	GST	CLK2	04-127	Full-length	1-499	NP_001281267.1	N-terminal GST	Insect (sf21)
CLK2	BTN	BTN-CLK2	04-427-20N	Full-length	1-499	NP_001281267.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	CLK3	04-128	Full-length	1-490	NP_003983.2	N-terminal GST	Insect (sf21)
CLK3	GST	CLK4	04-129	Full-length	1-481	NP_065717.1	N-terminal GST	Insect (sf21)
CLK4	BTN	BTN-CLK4	04-429-20N	Full-length	1-481	NP_065717.1	N-terminal DYKDDDDK	Insect (sf21)
	Tag-free	COT(MAP3K8)	07-301 (Tag-free)	Catalytic domain	30-397	NP_005195.2	Tag-free	Insect (sf21)
COT(MAP3K8)	GST	CRIK(CIT)	01-104	Catalytic domain	1-449	NP_009105.1	N-terminal GST	Insect (sf21)
CRIK(CIT)	GST	DAPK1	02-134	Catalytic domain	1-289	NP_004929.2	N-terminal GST	Insect (sf21)
DAPK1	GST	DAPK3	02-136	Full-length	1-454	NP_001339.1	N-terminal GST	Insect (sf21)
	BTN	BTN-DAPK3	02-436-20N	Full-length	1-454	NP_001339.1	N-terminal DYKDDDDK	Insect (sf21)
DAPK3	GST	DCAMKL1	02-139	Full-length	1-729	NP_004725.1	N-terminal GST	Insect (sf21)
DCAMKL1	GST	DCAMKL2	02-140	Truncated	1-691	NP_001035350.2	N-terminal GST	Insect (sf21)
DCAMKL2	GST	DLK(MAP3K12)	09-111	Catalytic domain	1-520	NP_006292.3	N-terminal GST	Insect (sf21)
	BTN	BTN-DLK(MAP3K12)	09-411-20N	Catalytic domain	1-520	NP_006292.3	N-terminal DYKDDDDK	Insect (sf21)
DLK(MAP3K12)	GST	DMPK1(DMPK)	01-105	Catalytic domain	1-544	NP_001075032.1	N-terminal GST	Insect (sf21)
DMPK1(DMPK)	GST	DMPK2(CDC42BPG)	01-106	Catalytic domain	1-473	NP_059995.2	N-terminal GST	Insect (sf21)
DMPK2(CDC42BPG)	GST	DRAK1(STK17A)	02-137	Full-length	1-414	NP_004751.2	N-terminal GST	Insect (sf21)
DRAK1(STK17A)	GST	DYRK1A	04-130	Full-length	1-763	NP_001387.2	N-terminal GST	Insect (sf21)
	BTN	BTN-DYRK1A	04-430-20N	Full-length	1-763	NP_001387.2	N-terminal DYKDDDDK	Insect (sf21)
DYRK1A	BTN							

Kinase Products

Serine / Threonine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
DYRK1B	GST	DYRK1B	04-131	Full-length	1-629	NP_004705.1	N-terminal GST	Insect (sf21)
DYRK2	GST	DYRK2	04-132	Full-length	1-528	NP_003574.1	N-terminal GST	Insect (sf21)
DYRK3	GST	DYRK3	04-133	Full-length	1-588	NP_003573.2	N-terminal GST	Insect (sf21)
DYRK4	DYKDDDDK	DYRK4	04-434	Full-length	1-520	NP_003836.1	N-terminal DYKDDDDK	Insect (sf21)
EEF2K	GST	EEF2K	10-113	Full-length	1-725	NP_037434.2	N-terminal GST	E. coli
Erk1(MAPK3)	GST	Erk1(MAPK3)	04-142	Full-length	1-379	NP_002737.2	N-terminal GST	E. coli
Erk2(MAPK1)	GST	Erk2(MAPK1)	04-143	Full-length	1-360	NP_002736.3	N-terminal GST	E. coli
Erk5(MAPK7)	GST	Erk5(MAPK7)	04-146	Catalytic domain	1-398	NP_002740.2	N-terminal GST	E. coli
Erk7(MAPK15)	GST	Erk7(MAPK15)	04-147	Full-length	1-544	NP_620590.2	N-terminal GST	Insect (sf21)
GCN2(EIF2AK4)	GST	GCN2(EIF2AK4)	05-153	Full-length	1-1649	NP_001013725	N-terminal GST	Insect (sf21)
	BTN	BTN-GCN2(EIF2AK4)	05-453-20N	Full-length	1-1649	NP_001013725	N-terminal DYKDDDDK	Insect (sf21)
GPRK4(GRK4)	Tag-free	GPRK4(GRK4)	01-313 (Tag-free)	Full-length	1-578	NP_892027.2	Tag-free	Insect (sf21)
GPRK6(GRK6)	Tag-free	GPRK6(GRK6)	01-315 (Tag-free)	Full-length	1-560	NP_001004105.2	Tag-free	Insect (sf21)
GPRK7(GRK7)	GST	GPRK7(GRK7)	01-116	Full-length	1-553	NP_631948.1	N-terminal GST	Insect (sf21)
GSK3α(GSK3A)	GST	GSK3α(GSK3A)	04-140	Full-length	1-483	NP_063937.2	N-terminal GST	Insect (sf21)
GSK3β(GSK3B)	GST	GSK3β(GSK3B)	04-141	Full-length	1-420	NP_001139628.1	N-terminal GST	Insect (sf21)
	BTN	BTN-GSK3β(GSK3B)	04-441-20N	Full-length	1-420	NP_001139628.1	N-terminal DYKDDDDK	Insect (sf21)
Haspin(GSG2)	GST	Haspin(GSG2)	05-111	Full-length	1-798	NP_114171.2	N-terminal GST	Insect (sf21)
HGK(MAP4K4)	GST	HGK(MAP4K4)	07-137	Catalytic domain	1-328	NP_004825.3	N-terminal GST	Insect (sf21)
HH498 (TNNI3K)	GST	HH498(TNNI3K)	09-112	Catalytic domain	1-727	NP_057062.1	N-terminal GST	Insect (sf21)
HIPK1	GST	HIPK1	04-135	Catalytic domain	158-555	NP_689909.2	N-terminal GST	Insect (sf21)
HIPK2	GST	HIPK2	04-136	Full-length	1-1198	NP_073577.3	N-terminal GST	Insect (sf21)
HIPK3	GST	HIPK3	04-137	Catalytic domain	161-562	NP_005725.3	N-terminal GST	Insect (sf21)
HIPK4	GST	HIPK4	04-138	Full-length	1-616	NP_653286.2	N-terminal GST	Insect (sf21)
HPK1(MAP4K1)	DYKDDDDK	HPK1(MAP4K1)	07-410	Catalytic domain	1-346	NP_009112.1	N-terminal DYKDDDDK	Insect (sf21)
HRI(EIF2AK1)	GST	HRI(EIF2AK1)	05-154	Full-length	1-630	NP_055228.2	N-terminal GST	E. coli
ICK	GST	ICK	04-156	Full-length	1-632	NP_057597.2	N-terminal GST	Insect (sf21)
IKKα(CHUK)	GST	IKKα(CHUK)	05-112	Full-length	1-745	NP_001269.3	N-terminal GST	Insect (sf21)
	BTN	BTN-IKKα(CHUK)	05-412-20N	Full-length	1-745	NP_001269.3	N-terminal DYKDDDDK	Insect (sf21)
IKKβ(IKBKB)	HIS	IKKβ(IKBKB)	05-084	Truncated	1-662	NP_001547.1	N-terminal His	Insect (sf21)
	BTN	BTN-IKKβ(IKBKB)	05-484-20N	Truncated	1-662	NP_001547.1	N-terminal DYKDDDDK	Insect (sf21)
IKKε(IKBKE)	GST	IKKε(IKBKE)	05-114	Full-length	1-716	NP_054721.1	N-terminal GST	Insect (sf21)
IRAK1	GST	IRAK1	09-101	Truncated	194-712	NP_001560.2	N-terminal GST	Insect (sf21)
IRAK4	GST	IRAK4	09-145	Full-length	1-460	NP_057207.2	N-terminal GST	Insect (sf21)
	BTN	BTN-IRAK4	09-445-20N	Full-length	1-460	NP_057207.2	N-terminal DYKDDDDK	Insect (sf21)
JNK1(MAPK8)	GST	JNK1(MAPK8)	04-163	Catalytic domain	2-364	NP_620634.1	N-terminal GST	E. coli
JNK2(MAPK9)	GST	JNK2(MAPK9)	04-164	Catalytic domain	1-364	NP_002743.3	N-terminal GST	E. coli
JNK3(MAPK10)	GST	JNK3(MAPK10)	04-150	Full-length	1-426	NP_620446.1	N-terminal GST	E. coli
LATS1/MOBKL1A	GST	LATS1/MOBKL1A	01-523	Catalytic domain	589-1130	NP_004681.1	N-terminal GST	Insect (sf21)
LATS2/MOBKL1A	GST	LATS2/MOBKL1A	01-524	Catalytic domain	553-1088	NP_055387.2	N-terminal GST	Insect (sf21)
	BTN	BTN-LATS2/MOBKL1A	01-424-20N	Catalytic domain	553-1088	NP_055387.2	N-terminal DYKDDDDK	Insect (sf21)
LIMK1	GST	LIMK1	09-105	Truncated	285-638	NP_002305.1	N-terminal GST	Insect (sf21)
	GST	LIMK2	09-106	Full-length	1-638	NP_005560.1	N-terminal GST	Insect (sf21)
LIMK2	BTN	BTN-LIMK2	09-406-20N	Full-length	1-638	NP_005560.1	N-terminal DYKDDDDK	Insect (sf21)
LKB1(STK11)/MO25α/STRADα	GST	LKB1(STK11)/MO25α/STRADα	02-119	Full-length	1-433	NP_000446.1	N-terminal GST	Insect (sf21)
LOK(STK10)	GST	LOK(STK10)	07-115	Full-length	1-968	NP_005981.3	N-terminal GST	Insect (sf21)
LRRK2	GST	LRRK2	09-110	Catalytic domain	960-2527	NP_940980.4	N-terminal GST	Insect (sf21)
LRRK2 [G2019S]	GST	LRRK2[G2019S]	09-146	Catalytic domain	960-2527	NP_940980.4	N-terminal GST	Insect (sf21)
LZK(MAP3K13)	GST	LZK(MAP3K13)	09-114	Full-length	1-966	NP_004712.1	N-terminal GST	Insect (sf21)
MAP2K1	GST	MAP2K1	07-141	Full-length	1-393	NP_002746.1	N-terminal GST	Insect (sf21)
MAP2K1 [F129L]	GST	MAP2K1[F129L]	07-151	Full-length	1-393	NP_002746.1	N-terminal GST	Insect (sf21)
MAP2K1 [P124L]	GST	MAP2K1[P124L]	07-150	Full-length	1-393	NP_002746.1	N-terminal GST	Insect (sf21)
MAP2K2	GST	MAP2K2	07-142	Full-length	1-400	NP_109587.1	N-terminal GST	Insect (sf21)
MAP2K3	GST	MAP2K3	07-143	Full-length	1-347	NP_659731.1	N-terminal GST	Insect (sf21)
MAP2K4	GST	MAP2K4	07-144	Full-length	1-399	NP_003001.1	N-terminal GST	Insect (sf21)
MAP2K5	GST	MAP2K5	07-145	Full-length	1-448	NP_660143.1	N-terminal GST	Insect (sf21)
MAP2K6	GST	MAP2K6	07-146	Full-length	1-334	NP_002749.2	N-terminal GST	Insect (sf21)
MAP2K7	GST	MAP2K7	07-148	Full-length	1-419	NP_660186.1	N-terminal GST	Insect (sf21)
MAP3K1	GST	MAP3K1	07-103	Catalytic domain	1327-1646	XP_042066.8	N-terminal GST	Insect (sf21)
MAP3K2	GST	MAP3K2	07-104	Catalytic domain	336-619	NP_006600.3	N-terminal GST	Insect (sf21)
MAP3K2(FL)	GST	MAP3K2(FL)	07-154	Full-length	1-619	NP_006600.3	N-terminal GST	Insect (sf21)
MAP3K3	GST	MAP3K3	07-105	Full-length	1-626	NP_002392.2	N-terminal GST	Insect (sf21)
MAP3K4	GST	MAP3K4	07-106	Catalytic domain	1312-1608	NP_005913.3	N-terminal GST	Insect (sf21)
MAP3K5	GST	MAP3K5	07-107	Catalytic domain	654-971	NP_005914.1	N-terminal GST	Insect (sf21)
	BTN	BTN-MAP3K5	07-407-20N	Catalytic domain	654-971	NP_005914.1	N-terminal DYKDDDDK	Insect (sf21)
MAP3K6	GST	MAP3K6	07-108	Catalytic domain	622-950	NP_004663.3	N-terminal GST	Insect (sf21)
MAP3K14	GST	MAP3K14	07-102	Catalytic domain	319-947	NP_003945.2	N-terminal GST	Insect (sf21)
MAP4K2	GST	MAP4K2	07-111	Full-length	1-820	NP_004570.2	N-terminal GST	Insect (sf21)
MAP4K5(KHS1)	GST	MAP4K5(KHS1)	07-113	Full-length	1-846	NP_006566.2	N-terminal GST	Insect (sf21)
MAPKAPK2	GST	MAPKAPK2	02-142	Full-length	1-400	NP_116584.2	N-terminal GST	Insect (sf21)
MAPKAPK3	GST	MAPKAPK3	02-143	Full-length	1-382	NP_004626.1	N-terminal GST	Insect (sf21)
MAPKAPK5	GST	MAPKAPK5	02-144	Full-length	1-471	NP_003659.2	N-terminal GST	Insect (sf21)
MARK1	GST	MARK1	02-120	Full-length	1-795	NP_061120.3	N-terminal GST	Insect (sf21)
MARK2	GST	MARK2	02-121	Full-length	1-745	NP_059672.2	N-terminal GST	Insect (sf21)
MARK3	GST	MARK3	02-122	Full-length	1-729	NP_002367.5	N-terminal GST	Insect (sf21)
MARK4	GST	MARK4	02-123	Full-length	1-688	NP_113605.2	N-terminal GST	Insect (sf21)
MELK	GST	MELK	02-124	Truncated	1-493	NP_055606.1	N-terminal GST	E. coli
MINK(MINK1)	GST	MINK(MINK1)	07-139	Catalytic domain	1-314	NP_056531.1	N-terminal GST	Insect (sf21)
MLK1(MAP3K9)	GST	MLK1(MAP3K9)	09-115	Catalytic domain	110-422	NP_149132.2	N-terminal GST	Insect (sf21)
	BTN	BTN-MLK1(MAP3K9)	09-415-20N	Catalytic domain	110-422	NP_149132.2	N-terminal DYKDDDDK	Insect (sf21)

Kinase Products

Serine / Threonine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
MLK2(MAP3K10)	GST	MLK2(MAP3K10)	09-116	Catalytic domain	75-462	NP_002437.2	N-terminal GST	Insect (sf21)
	BTN	BTN-MLK2(MAP3K10)	09-416-20N	Catalytic domain	75-462	NP_002437.2	N-terminal DYKDDDDK	Insect (sf21)
MLK3(MAP3K11)	GST	MLK3(MAP3K11)	09-117	Catalytic domain	99-398	NP_002410.1	N-terminal GST	Insect (sf21)
	BTN	BTN-MLK3(MAP3K11)	09-417-20N	Catalytic domain	99-398	NP_002410.1	N-terminal DYKDDDDK	Insect (sf21)
MLK4	GST	MLK4	09-118	Catalytic domain	114-420	NP_115811.2	N-terminal GST	Insect (sf21)
MNK1(MKNK1)	GST	MNK1(MKNK1)	02-145	Full-length	1-424	NP_001129025.1	N-terminal GST	Insect (sf21)
	BTN	BTN-MNK1(MKNK1)	02-445-20N	Full-length	1-424	NP_001129025.1	N-terminal DYKDDDDK	Insect (sf21)
MNK2(MKNK2)	GST	MNK2(MKNK2)	02-146	Full-length	1-465	NP_951009.1	N-terminal GST	Insect (sf21)
MOS	GST	MOS	05-118	Full-length	1-346	NP_005363.1	N-terminal GST	Insect (sf21)
MRCKα(CDC42BPA)	GST	MRCKα(CDC42BPA)	01-107	Truncated	1-574	NP_003598.2	N-terminal GST	Insect (sf21)
MRCKβ(CDC42BPB)	GST	MRCKβ(CDC42BPB)	01-108	Truncated	1-473	NP_006026.3	N-terminal GST	Insect (sf21)
MSK1(RPS6KA5)	GST	MSK1(RPS6KA5)	01-147	Full-length	1-802	NP_004746.2	N-terminal GST	Insect (sf21)
MSK2(RPS6KA4)	GST	MSK2(RPS6KA4)	01-148	Full-length	1-772	NP_003933.1	N-terminal GST	Insect (sf21)
MSSK1(STK23)	GST	MSSK1(STK23)	04-159	Full-length	1-567	NP_055185.2	N-terminal GST	Insect (sf21)
MST1(STK4)	GST	MST1(STK4)	07-116	Full-length	1-487	NP_006273.1	N-terminal GST	Insect (sf21)
MST2(STK3)	GST	MST2(STK3)	07-117	Full-length	1-491	NP_006272.2	N-terminal GST	Insect (sf21)
	BTN	BTN-MST2(STK3)	07-417-20N	Full-length	1-491	NP_006272.2	N-terminal DYKDDDDK	Insect (sf21)
MST3(STK24)	GST	MST3(STK24)	07-118	Full-length	1-431	NP_001027467.2	N-terminal GST	Insect (sf21)
MST4	GST	MST4	07-119	Full-length	1-416	NP_057626.2	N-terminal GST	Insect (sf21)
MYLK4	GST	MYLK4	02-161	Full-length	1-388	NP_001012418.2	N-terminal GST	Insect (sf21)
	BTN	BTN-MYLK4	02-461-20N	Full-length	1-388	NP_001012418.2	N-terminal DYKDDDDK	Insect (sf21)
MYT1(PKMYT1)	GST	MYT1(PKMYT1)	05-176	Full-length	1-499	NP_004194.3	N-terminal GST	Insect (sf21)
NDRI(STK38)	GST	NDRI(STK38)	01-125	Full-length	1-465	NP_009202.1	N-terminal GST	Insect (sf21)
NDR2(STK38L)	GST	NDR2(STK38L)	01-126	Full-length	1-464	NP_055815.1	N-terminal GST	Insect (sf21)
NEK1	GST	NEK1	05-123	Catalytic domain	1-505	NP_036356.1	N-terminal GST	Insect (sf21)
NEK2	Tag-free	NEK2	05-226 (Tag-free)	Full-length	1-445	NP_002488.1	Tag-free	Insect (sf21)
NEK3	GST	NEK3	05-127	Full-length	1-506	NP_002489.1	N-terminal GST	Insect (sf21)
NEK4	GST	NEK4	05-128	Full-length	1-841	NP_003148.2	N-terminal GST	Insect (sf21)
NEK6	GST	NEK6	05-130	Full-length	1-313	NP_055212.2	N-terminal GST	Insect (sf21)
NEK7	GST	NEK7	05-131	Full-length	1-302	NP_598001.1	N-terminal GST	Insect (sf21)
NEK9	GST	NEK9	05-133	Truncated	1-346, 733-979	NP_149107.4	N-terminal GST	Insect (sf21)
	BTN	BTN-NEK9	05-433-20N	Truncated	1-346, 733-979	NP_149107.4	N-terminal DYKDDDDK	Insect (sf21)
NEK11	GST	NEK11	05-125	Full-length	1-645	NP_079076.3	N-terminal GST	Insect (sf21)
NIM1K(MGC42105)	GST	NIM1K(MGC42105)	02-175	Full-length	1-436	NP_699192.1	N-terminal GST	Insect (sf21)
NLK	GST	NLK	04-151	Truncated	133-527	NP_057315.3	N-terminal GST	Insect (sf21)
NuaK1	GST	NuaK1	02-126	Full-length	1-661	NP_055655.1	N-terminal GST	Insect (sf21)
NuaK2	GST	NuaK2	02-127	Full-length	1-628	NP_112214.3	N-terminal GST	Insect (sf21)
	BTN	BTN-NuaK2	02-427-20N	Full-length	1-628	NP_112214.3	N-terminal DYKDDDDK	Insect (sf21)
OSR1(OXSR1)	GST	OSR1(OXSR1)	07-122	Full-length	1-527	NP_005100.1	N-terminal GST	Insect (sf21)
p38α(MAPK14)	GST	p38α(MAPK14)	04-152	Truncated	9-352	NP_620581.1	N-terminal GST	E. coli
	BTN	BTN-p38α(MAPK14)	04-452-20N	Truncated	9-352	NP_620581.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-p38α(MAPK14)[inactive]	04-452-10-20N	Truncated	9-352	NP_620581.1	N-terminal DYKDDDDK	Insect (sf21)
p38β(MAPK11)	GST	p38β(MAPK11)	04-153	Full-length	1-364	NP_002742.3	N-terminal GST	E. coli
p38γ(MAPK12)	GST	p38γ(MAPK12)	04-155	Full-length	1-367	NP_002960.2	N-terminal GST	E. coli
p38δ(MAPK13)	GST	p38δ(MAPK13)	04-154	Full-length	1-365	NP_002745.1	N-terminal GST	E. coli
p70S6K(RPS6KB1)	GST	p70S6K(RPS6KB1)	01-156	Catalytic domain	1-421	NP_003152.1	N-terminal GST	Insect (sf21)
p70S6Kβ(RPS6KB2)	GST	p70S6Kβ(RPS6KB2)	01-155	Full-length	1-482	NP_003943.2	N-terminal GST	Insect (sf21)
PAK1	GST	PAK1	07-123	Full-length	1-545	NP_002567.3	N-terminal GST	Insect (sf21)
PAK2	GST	PAK2	07-124	Full-length	1-524	NP_002568.2	N-terminal GST	Insect (sf21)
PAK3	GST	PAK3	07-125	Full-length	1-544	NP_002569.1	N-terminal GST	Insect (sf21)
PAK4	GST	PAK4	07-126	Full-length	1-591	NP_005875.1	N-terminal GST	Insect (sf21)
	BTN	BTN-PAK4	07-426-20N	Full-length	1-591	NP_005875.1	N-terminal DYKDDDDK	Insect (sf21)
PAK5(PAK7)	GST	PAK5(PAK7)	07-127	Catalytic domain	425-719	NP_065074.1	N-terminal GST	Insect (sf21)
PAK6	GST	PAK6	07-128	Full-length	1-681	NP_064553.1	N-terminal GST	Insect (sf21)
	BTN	BTN-PAK6	07-428-20N	Full-length	1-681	NP_064553.1	N-terminal DYKDDDDK	Insect (sf21)
PASK	GST	PASK	02-128	Catalytic domain	949-1323	NP_055963.2	N-terminal GST	Insect (sf21)
PBK	GST	PBK	05-168	Full-length	1-322	NP_060962.2	N-terminal GST	Insect (sf21)
PCTAIRE1(CDK16)/CycY	GST	PCTAIRE1(CDK16)/CycY	04-116	Catalytic domain	107-496	NP_006192.1	N-terminal GST	Insect (sf21)
PDHK1(PDK1)	GST	PDHK1(PDK1)	10-123	Full-length	1-436	NP_002601.1	N-terminal GST	Insect (sf21)
PDHK2(PDK2)	GST	PDHK2(PDK2)	10-140	Full-length	1-407	NP_002602.2	N-terminal GST	Insect (sf21)
PDHK3(PDK3)	GST	PDHK3(PDK3)	10-124	Full-length	1-406	NP_005382.1	N-terminal GST	Insect (sf21)
PDHK4(PDK4)	GST	PDHK4(PDK4)	10-125	Full-length	1-411	NP_002603.1	N-terminal GST	E. coli
PDK1(PDPK1)	GST	PDK1(PDPK1)	01-132	Full-length	1-556	NP_002604.1	N-terminal GST	Insect (sf21)
PEK(EIF2AK3)	GST	PEK(EIF2AK3)	05-155	Cytoplasmic domain	536-1116	NP_004827.4	N-terminal GST	E. coli
PGK(PRKG1)	GST	PGK(PRKG1)	01-142	Full-length	1-686	NP_006249.1	N-terminal GST	Insect (sf21)
PHKG1	GST	PHKG1	02-152	Full-length	1-387	NP_006204.1	N-terminal GST	Insect (sf21)
PHKG2	GST	PHKG2	02-153	Full-length	1-406	NP_000285.1	N-terminal GST	Insect (sf21)
PIM1	HIS	PIM1	02-054	Full-length	1-313	NP_002639.1	N-terminal His	Insect (sf21)
PIM2	GST	PIM2	02-155	Full-length	1-311	NP_006866.2	N-terminal GST	Insect (sf21)
PIM3	GST	PIM3	02-156	Full-length	1-326	NP_001001852.2	N-terminal GST	Insect (sf21)
PKAcα(PRKACA)	GST	PKAcα(PRKACA)	01-127	Full-length	1-351	NP_002721.1	N-terminal GST	Insect (sf21)
	BTN	BTN-PKAcα(PRKACA)	01-427-20N	Full-length	1-351	NP_002721.1	N-terminal DYKDDDDK	Insect (sf21)
PKACβ(PRKACB)	GST	PKACβ(PRKACB)	01-128	Full-length	1-351	NP_002722.1	N-terminal GST	Insect (sf21)
PKACγ(PRKACG)	GST	PKACγ(PRKACG)	01-129	Full-length	1-351	NP_002723.2	N-terminal GST	Insect (sf21)
PKCα(PRKCA)	GST	PKCα(PRKCA)	01-133	Full-length	1-672	NP_002728.2	N-terminal GST	Insect (sf21)
	BTN	BTN-PKCα(PRKCA)	01-433-20N	Full-length	1-672	NP_002728.2	N-terminal DYKDDDDK	Insect (sf21)
PKCβ1(PRKCB1)	GST	PKCβ1(PRKCB1)	01-134	Full-length	1-671	NP_997700.1	N-terminal GST	Insect (sf21)
PKCβ2(PRKCB2)	GST	PKCβ2(PRKCB2)	01-165	Full-length	1-673	NP_002729.2	N-terminal GST	Insect (sf21)
PKCγ(PRKCG)	GST	PKCγ(PRKCG)	01-137	Full-length	1-697	NP_002730.1	N-terminal GST	Insect (sf21)
	BTN	BTN-PKCγ(PRKCG)	01-437-20N	Full-length	1-697	NP_002730.1	N-terminal DYKDDDDK	Insect (sf21)

Kinase Products

Serine / Threonine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	Carna Product Name	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
PKCδ(PRKCD)	GST	PKCδ(PRKCD)	01-135	Full-length	1-676	NP_006245.2	N-terminal GST	Insect (sf21)
PKCε(PRKCE)	GST	PKCε(PRKCE)	01-136	Full-length	1-737	NP_005391.1	N-terminal GST	Insect (sf21)
PKCζ(PRK CZ)	GST	PKCζ(PRK CZ)	01-141	Full-length	1-592	NP_002735.3	N-terminal GST	Insect (sf21)
PKCη(PRKCH)	GST	PKCη(PRKCH)	01-138	Full-length	1-683	NP_006246.2	N-terminal GST	Insect (sf21)
	BTN	BTN-PKCη(PRKCH)	01-438-20N	Full-length	1-683	NP_006246	N-terminal DYKDDDDK	Insect (sf21)
PKCθ(PRK CQ)	GST	PKCθ(PRK CQ)	01-140	Full-length	1-706	NP_006248.1	N-terminal GST	Insect (sf21)
	BTN	BTN-PKCθ(PRK CQ)	01-440-20N	Full-length	1-706	NP_006248.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	BTN-PKCθ(PRK CQ)[non-activated]	01-440-23N	Full-length	1-706	NP_006248.1	N-terminal DYKDDDDK	Insect (sf21)
	GST	PKCθ(PRKCI)	01-139	Truncated	10-596	NP_002731.4	N-terminal GST	Insect (sf21)
PKN1	GST	PKN1	01-144	Full-length	1-942	NP_002732.3	N-terminal GST	Insect (sf21)
	BTN	BTN-PKN1	01-444-20N	Full-length	1-942	NP_002732.3	N-terminal DYKDDDDK	Insect (sf21)
PKN2	GST	PKN2	01-145	Full-length	1-984	NP_006247.1	N-terminal GST	Insect (sf21)
PKN3	GST	PKN3	01-146	Full-length	1-889	NP_037487.2	N-terminal GST	Insect (sf21)
PKR(EIF2AK2)	GST	PKR(EIF2AK2)	05-156	Catalytic domain	252-551	NP_002750.1	N-terminal GST	Insect (sf21)
PLK1	GST	PLK1	05-157	Full-length	1-603	NP_005021.2	N-terminal GST	Insect (sf21)
	BTN	BTN-PLK1	05-457-20N	Full-length	1-603	NP_005021.2	N-terminal DYKDDDDK	Insect (sf21)
PLK2	GST	PLK2	05-158	Full-length	1-685	NP_006613.2	N-terminal GST	Insect (sf21)
PLK3	GST	PLK3	05-159	Catalytic domain	58-340	NP_004064.2	N-terminal GST	Insect (sf21)
PRKD1(PKD1)	GST	PRKD1(PKD1)	02-157	Full-length	1-912	NP_002733.2	N-terminal GST	Insect (sf21)
PRKD2(PKD2)	GST	PRKD2(PKD2)	02-158	Full-length	1-878	NP_057541.2	N-terminal GST	Insect (sf21)
	BTN	BTN-PRKD2(PKD2)	02-458-20N	Full-length	1-878	NP_057541.2	N-terminal DYKDDDDK	Insect (sf21)
PRKD3(PKD3)	GST	PRKD3(PKD3)	02-159	Full-length	1-890	NP_005804.1	N-terminal GST	Insect (sf21)
PRKX	GST	PRKX	01-130	Full-length	1-358	NP_005035.1	N-terminal GST	Insect (sf21)
QIK(SNF1LK2)	GST	QIK(SNF1LK2)	02-129	Full-length	1-926	NP_056006.1	N-terminal GST	Insect (sf21)
RAF1	GST	RAF1	09-125	Catalytic domain	306-648	NP_002871.1	N-terminal GST	Insect (sf21)
	BTN	BTN-RAF1	09-425-20N	Catalytic domain	306-648	NP_002871.1	N-terminal DYKDDDDK	Insect (sf21)
RIPK2	GST	RIPK2	09-128	Catalytic domain	1-305	NP_003812.1	N-terminal GST	Insect (sf21)
ROCK1	GST	ROCK1	01-109	Catalytic domain	1-477	NP_005397.1	N-terminal GST	Insect (sf21)
	BTN	BTN-ROCK1	01-409-20N	Catalytic domain	1-477	NP_005397.1	N-terminal DYKDDDDK	Insect (sf21)
ROCK2	GST	ROCK2	01-110	Catalytic domain	1-553	NP_004841.2	N-terminal GST	Insect (sf21)
	BTN	BTN-ROCK2	01-410-20N	Catalytic domain	1-553	NP_004841.2	N-terminal DYKDDDDK	Insect (sf21)
RSK1(RPS6KA1)	GST	RSK1(RPS6KA1)	01-149	Full-length	1-735	NP_002944.2	N-terminal GST	Insect (sf21)
RSK2(RPS6KA3)	GST	RSK2(RPS6KA3)	01-150	Full-length	1-740	NP_004577.1	N-terminal GST	Insect (sf21)
RSK3(RPS6KA2)	GST	RSK3(RPS6KA2)	01-151	Full-length	1-733	NP_066958.2	N-terminal GST	Insect (sf21)
RSK4(RPS6KA6)	GST	RSK4(RPS6KA6)	01-152	Full-length	1-745	NP_055311.1	N-terminal GST	Insect (sf21)
SGK	GST	SGK	01-158	Truncated	61-431	NP_005618.2	N-terminal GST	Insect (sf21)
SGK2	GST	SGK2	01-159	Full-length	1-367	NP_733794.1	N-terminal GST	Insect (sf21)
SGK3(SGKL)	GST	SGK3(SGKL)	01-160	Truncated	119-496	NP_037389.4	N-terminal GST	Insect (sf21)
SIK(SNF1LK)	GST	SIK(SNF1LK)	02-131	Full-length	1-783	NP_775490.2	N-terminal GST	Insect (sf21)
SIK3(QSK)	GST	SIK3(QSK)	02-130	Full-length	1-1369	NP_001353615.1	N-terminal GST	Insect (sf21)
SIK3(QSK)(59-365aa)	GST	SIK3(QSK)(59-365aa)	02-133	Catalytic domain	59-365	NP_001353615.1	N-terminal GST	Insect (sf21)
skMLCK(MYLK2)	GST	skMLCK(MYLK2)	02-150	Full-length	1-596	NP_149109.1	N-terminal GST	Insect (sf21)
SLK	GST	SLK	07-129	Full-length	1-1152	NP_055535.1	N-terminal GST	Insect (sf21)
smMLCK(MYLK)	GST	smMLCK(MYLK)	02-151	Catalytic domain	1428-1771	NP_444253.3	N-terminal GST	Insect (sf21)
SRPK1	GST	SRPK1	04-160	Full-length	1-655	NP_003128.3	N-terminal GST	E. coli
	BTN	BTN-SRPK1	04-460-20N	Full-length	1-655	NP_003128.3	N-terminal DYKDDDDK	Insect (sf21)
SRPK2	GST	SRPK2	04-161	Full-length	1-688	NP_872633.1	N-terminal GST	Insect (sf21)
SSTK(TSSK6)	GST	SSTK(TSSK6)	02-163	Full-length	1-273	NP_114426.1	N-terminal GST	Insect (sf21)
STK33	Tag-free	STK33	02-301 (Tag-free)	Full-length	1-514	NP_112168.1	Tag-free	Insect (sf21)
STLK3(STK39)	GST	STLK3(STK39)	07-130	Full-length	1-545	NP_037365.2	N-terminal GST	Insect (sf21)
TAK1-TAB1(MAP3K7)	HIS	TAK1-TAB1(MAP3K7)	09-019	Fused	1-303	NP_663304.1	N-terminal His	Insect (sf21)
	BTN	BTN-TAK1-TAB1(MAP3K7)	09-419-21C	Fused	1-303	NP_663304.1	C-terminal DYKDDDDK	Insect (sf21)
TAOK2	GST	TAOK2	07-133	Catalytic domain	1-319	NP_004774.1	N-terminal GST	Insect (sf21)
TAOK3	GST	TAOK3	07-135	Full-length	1-898	NP_057365.3	N-terminal GST	Insect (sf21)
TBK1	GST	TBK1	05-115	Full-length	1-729	NP_037386.1	N-terminal GST	Insect (sf21)
	BTN	BTN-TBK1	05-415-20N	Full-length	1-729	NP_037386.1	N-terminal DYKDDDDK	Insect (sf21)
TESK1	GST	TESK1	09-107	Full-length	1-626	NP_006276.2	N-terminal GST	Insect (sf21)
TGFβR1(TGFBR1)	GST	TGFβR1(TGFBR1)	09-141	Catalytic domain	200-503	NP_004603.1	N-terminal GST	Insect (sf21)
	BTN	BTN-TGFβR1(TGFBR1)	09-441-20N	Catalytic domain	200-503	NP_004603.1	N-terminal DYKDDDDK	Insect (sf21)
TGFβR2(TGFBR2)	GST	TGFβR2(TGFBR2)	09-142	Catalytic domain	194-567	NP_003233.4	N-terminal GST	Insect (sf21)
TLK1	GST	TLK1	05-166	Full-length	1-766	NP_036422.3	N-terminal GST	Insect (sf21)
TLK2	GST	TLK2	05-167	Full-length	1-750	NP_006843.2	N-terminal GST	Insect (sf21)
TNIK	GST	TNIK	07-138	Catalytic domain	1-314	NP_055843.1	N-terminal GST	Insect (sf21)
	BTN	BTN-TNIK	07-438-20N	Catalytic domain	1-314	NP_055843.1	N-terminal DYKDDDDK	Insect (sf21)
TSSK1B(TSSK1)	Tag-free	TSSK1B(TSSK1)	02-364 (Tag-free)	Full-length	1-367	NP_114417.1	Tag-free	Insect (sf21)
	BTN	BTN-TSSK1B(TSSK1)	02-464-20N	Full-length	1-367	NP_114417.1	N-terminal DYKDDDDK	Insect (sf21)
TSSK2	GST	TSSK2	02-165	Full-length	1-358	NP_443732.3	N-terminal GST	Insect (sf21)
	BTN	BTN-TSSK2	02-465-20N	Full-length	1-358	NP_443732.3	N-terminal DYKDDDDK	Insect (sf21)
TSSK3	GST	TSSK3	02-166	Full-length	2-268	NP_443073.1	N-terminal GST	Insect (sf21)
TTBK1	GST	TTBK1	03-108	Catalytic domain	1-343	NP_115927.1	N-terminal GST	Insect (sf21)
TTBK2	GST	TTBK2	03-109	Catalytic domain	1-331	NP_775771.3	N-terminal GST	Insect (sf21)
TTK	GST	TTK	05-169	Full-length	1-857	NP_003309.2	N-terminal GST	Insect (sf21)

Kinase Products

Serine / Threonine Kinases

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	CarnaProductName	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
ULK3	GST	ULK3	05-173	Full-length	1-472	NP_001092906.3	N-terminal GST	Insect (sf21)
VRK1	GST	VRK1	03-110	Full-length	1-396	NP_003375.1	N-terminal GST	E. coli
VRK2	GST	VRK2	03-111	Catalytic domain	1-483	NP_006287.2	N-terminal GST	Insect (sf21)
WEE1	GST	WEE1	05-177	Catalytic domain	215-646	NP_003381.1	N-terminal GST	Insect (sf21)
	BTN	BTN-WEE1	05-477-20N	Catalytic domain	215-646	NP_003381.1	N-terminal DYKDDDDK	Insect (sf21)
WNK1	GST	WNK1	05-179	Catalytic domain	1-491	NP_061852.3	N-terminal GST	Insect (sf21)
	BTN	BTN-WNK1	05-479-20N	Catalytic domain	1-491	NP_061852.3	N-terminal DYKDDDDK	Insect (sf21)
WNK2	GST	WNK2	05-180	Catalytic domain	166-489	NP_006639.3	N-terminal GST	Insect (sf21)
WNK3	GST	WNK3	05-181	Catalytic domain	1-434	NP_065973.2	N-terminal GST	Insect (sf21)
WNK4	GST	WNK4	05-182	Catalytic domain	1-444	NP_115763.2	N-terminal GST	Insect (sf21)
	BTN	BTN-WNK4	05-482-20N	Catalytic domain	1-444	NP_115763.2	N-terminal DYKDDDDK	Insect (sf21)
YSK1(STK25)	GST	YSK1(STK25)	07-136	Full-length	1-426	NP_006365.2	N-terminal GST	Insect (sf21)
ZAK	GST	ZAK	09-120	Full-length	1-455	NP_598407.1	N-terminal GST	Insect (sf21)

Lipid Kinases

Protein Kinase	Tag	CarnaProductName	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
DGKα(DGKA)	GST	DGKα(DGKA)	12-101	Full-length	1-735	NP_958852.1	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKα(DGKA)	12-401-20N	Full-length	1-735	NP_958852.1	N-terminal DYKDDDDK	Insect (sf21)
DGKα(DGKA)(196-735aa)	BTN	BTN-DGKα(DGKA)(196-735aa)	12-411-20N	Truncated	196-735	NP_958852.1	N-terminal DYKDDDDK	Insect (sf21)
DGKβ(DGKB)	GST	DGKβ(DGKB)	12-102	Full-length	1-803	NP_001337638.1	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKβ(DGKB)	12-402-20N	Full-length	1-803	NP_001337638.1	N-terminal DYKDDDDK	Insect (sf21)
	BTN	mouse BTN-DGKβ(DGKB)	12M-402-20N	Full-length	1-802	NP_001348615.1	N-terminal DYKDDDDK	Insect (sf21)
DGKβ(DGKB)(235-803aa)	BTN	BTN-DGKβ(DGKB)(235-803aa)	12-412-20N	Truncated	235-803	NP_001337638.1	N-terminal DYKDDDDK	Insect (sf21)
DGKγ(DGKG)	GST	DGKγ(DGKG)	12-103	Full-length	1-791	NP_001337.2	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKγ(DGKG)	12-403-20N	Full-length	1-791	NP_001337.2	N-terminal DYKDDDDK	Insect (sf21)
	BTN	mouse BTN-DGKγ(DGKG)	12M-403-20N	Full-length	1-788	NP_619591.1	N-terminal DYKDDDDK	Insect (sf21)
DGKγ(DGKG)(261-791aa)	BTN	BTN-DGKγ(DGKG)(261-791aa)	12-413-20N	Truncated	261-791	NP_001337.2	N-terminal DYKDDDDK	Insect (sf21)
DGKδ(DGKD)	GST	DGKδ(DGKD)	12-104	Catalytic domain	1-1141	NP_690618.2	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKδ(DGKD)	12-404-20N	Catalytic domain	1-1141	NP_690618.2	N-terminal DYKDDDDK	Insect (sf21)
DGKε(DGKE)	GST	DGKε(DGKE)	12-115	Truncated protein	48-567	NP_003638.1	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKε(DGKE)	12-415-20N	Truncated protein	48-567	NP_003638.1	N-terminal DYKDDDDK	Insect (sf21)
DGKη(DGKH)	GST	DGKη(DGKH)	12-106	Catalytic domain	1-1147	NP_821077.1	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKη(DGKH)	12-406-20N	Catalytic domain	1-1147	NP_821077.1	N-terminal DYKDDDDK	Insect (sf21)
DGKι(DGKI)	GST	DGKι(DGKI)	12-107	Full-length	1-1065	NP_004708.1	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKι(DGKI)	12-407-20N	Full-length	1-1065	NP_004708.1	N-terminal DYKDDDDK	Insect (sf21)
DGKκ(DGKK)	GST	DGKκ(DGKK)	12-108	Catalytic domain	216-1271	NP_001013764.1	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKκ(DGKK)	12-408-20N	Catalytic domain	216-1271	NP_001013764.1	N-terminal DYKDDDDK	Insect (sf21)
DGKθ(DGKQ)	GST	DGKθ(DGKQ)	12-109	Full-length	1-942	NP_001338.2	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKθ(DGKQ)	12-409-20N	Full-length	1-942	NP_001338.2	N-terminal DYKDDDDK	Insect (sf21)
DGKζ(DGKZ)	GST	DGKζ(DGKZ)	12-110	Full-length	1-929	NP_003637.2	N-terminal GST	Insect (sf21)
	BTN	BTN-DGKζ(DGKZ)	12-410-20N	Full-length	1-929	NP_003637.2	N-terminal DYKDDDDK	Insect (sf21)
MTOR(FRAP)/MLST8	DYKDDDDK	MTOR(FRAP)/MLST8	11-431	Catalytic domain	1362-2549(MTOR), 1-326(MLST8)	NP_004949.1(MTOR), NP_071767.3(MLST8)	N-terminal DYKDDDDK	Insect (sf21)
PI4KB	GST	PI4KB	11-131	Full-length	1-828	NP_002642.1	N-terminal GST	Insect (sf21)
PIK3C3	GST	PIK3C3	11-122	Full-length	1-887	NP_002638.2	N-terminal GST	Insect (sf21)
PIK3CA/PIK3R1	GST	PIK3CA/PIK3R1	11-101	Full-length	1-1068	NP_006209.2	N-terminal GST	Insect (sf21)
	BTN	BTN-PIK3CA/PIK3R1	11-401-20N	Full-length	1-1068	NP_006209.2	N-terminal DYKDDDDK	Insect (sf21)
PIK3CA [E542K]/PIK3R1	BTN	BTN-PIK3CA[E542K]/PIK3R1	11-413-20N	Full-length	1-1068	NP_006209.2	N-terminal DYKDDDDK	Insect (sf21)
PIK3CA [E545K]/PIK3R1	BTN	BTN-PIK3CA[E545K]/PIK3R1	11-414-20N	Full-length	1-1068	NP_006209.2	N-terminal DYKDDDDK	Insect (sf21)
PIK3CA [H1047R]/PIK3R1	BTN	BTN-PIK3CA[H1047R]/PIK3R1	11-415-20N	Full-length	1-1068	NP_006209.2	N-terminal DYKDDDDK	Insect (sf21)
PIK3CA [P539R]/PIK3R1	BTN	BTN-PIK3CA[P539R]/PIK3R1	11-412-20N	Full-length	1-1068	NP_006209.2	N-terminal DYKDDDDK	Insect (sf21)
PIK3CA [R88Q]/PIK3R1	BTN	BTN-PIK3CA[R88Q]/PIK3R1	11-411-20N	Full-length	1-1068	NP_006209.2	N-terminal DYKDDDDK	Insect (sf21)
PIK3CB/PIK3R1	GST	PIK3CB/PIK3R1	11-102	Full-length	1-1070	NP_006210.1	N-terminal GST	Insect (sf21)
	BTN	BTN-PIK3CB/PIK3R1	11-402-20N	Full-length	1-1070	NP_006210.1	N-terminal DYKDDDDK	Insect (sf21)
PIK3CD/PIK3R1	GST	PIK3CD/PIK3R1	11-103	Full-length	1-1044	NP_005017.3	N-terminal GST	Insect (sf21)
	BTN	BTN-PIK3CD/PIK3R1	11-403-20N	Full-length	1-1044	NP_005017.3	N-terminal DYKDDDDK	Insect (sf21)
PIK3CG	BTN	BTN-PIK3CG	11-404-20N	Full-length	1-1102	NP_002640.2	N-terminal DYKDDDDK	Insect (sf21)
PIKFYVE(PIPSK3)	GST	PIKFYVE(PIPSK3)	11-118	Full-length	1-2098	NP_055855.2	N-terminal GST	Insect (sf21)
	BTN	BTN-PIKFYVE(PIPSK3)	11-418-20N	Full-length	1-2098	NP_055855.2	N-terminal DYKDDDDK	Insect (sf21)
PIP4K2A	GST	PIP4K2A	11-115	Full-length	1-406	NP_005019	N-terminal GST	Insect (sf21)
PIP4K2B	GST	PIP4K2B	11-116	Full-length	1-416	NP_003550	N-terminal GST	Insect (sf21)
PIP4K2C	GST	PIP4K2C	11-117	Full-length	1-421	NP_079055	N-terminal GST	Insect (sf21)
PIP5K1A	GST	PIP5K1A	11-111	Full-length	1-549	NP_003548.1	N-terminal GST	Insect (sf21)
PIP5K1B	GST	PIP5K1B	11-112	Full-length	1-540	NP_003549.1	N-terminal GST	Insect (sf21)
PIP5K1C	GST	PIP5K1C	11-113	Full-length	1-668	NP_036530	N-terminal GST	Insect (sf21)
PIP5KL1	GST	PIP5KL1	11-114	Full-length	1-394	NP_001128691.1	N-terminal GST	Insect (sf21)
SPHK1	GST	SPHK1	11-105	Full-length	1-384	NP_001136074.1	N-terminal GST	Insect (sf21)
	BTN	BTN-SPHK1	11-405-20N	Full-length	1-384	NP_001136074.1	N-terminal DYKDDDDK	Insect (sf21)
SPHK2	GST	SPHK2	11-106	Full-length	1-618	NP_001191089.1	N-terminal GST	Insect (sf21)
	BTN	BTN-SPHK2	11-406-20N	Full-length	1-618	NP_001191089.1	N-terminal DYKDDDDK	Insect (sf21)

Kinase Products

Inactive Kinases**

For details of products, please see "PRODUCTS AND SERVICES" on page 1~3.
Note: Please contact us for availability or further information. Information may be changed without notice.

Protein Kinase	Tag	CarnaProductName	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
Erk2(MAPK1) [inactive]	GST	Erk2(MAPK1)[inactive]	04-143-10	Full-length	1-360	NP_002736.3	N-terminal GST	E. coli
Erk5(MAPK7) [inactive]	GST	Erk5(MAPK7)[inactive]	04-146-10	Catalytic domain	1-398	NP_002740.2	N-terminal GST	E. coli
JNK1(MAPK8) [inactive]	GST	JNK1(MAPK8)[inactive]	04-163-10	Catalytic domain	2-364	NP_620634.1	N-terminal GST	E. coli
JNK2(MAPK9) [inactive]	GST	JNK2(MAPK9)[inactive]	04-164-10	Catalytic domain	1-364	NP_002743.3	N-terminal GST	E. coli
MAP2K1 [inactive]	GST	MAP2K1[inactive]	07-141-10	Full-length	1-393	NP_002746.1	N-terminal GST	E. coli
	BTN	BTN-MAP2K1[inactive]	07-441-10-20N	Full-length	1-393	NP_002746.1	N-terminal DYKDDDDK	Insect (sf21)
MAP2K6 [inactive]	GST	MAP2K6[inactive]	07-146-10	Full-length	1-334	NP_002749.2	N-terminal GST	E. coli
MAP2K7 [inactive]	GST	MAP2K7[inactive]	07-147-10	Full-length	1-419	NP_660186.1	N-terminal GST	E. coli
p38α(MAPK14) [inactive]	GST	p38α(MAPK14)[inactive]	04-152-10	Truncated	9-352	NP_620581.1	N-terminal GST	E. coli

Inactive Mutant Kinases**

Protein Kinase	Tag	CarnaProductName	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
Erk2(MAPK1) [inactive mutant]	GST	Erk2(MAPK1)[inactivemutant]	04-143-11	Full-length	1-360	NP_002736.3	N-terminal GST	E. coli
Erk5(MAPK7) [inactive mutant]	GST	Erk5(MAPK7)[inactivemutant]	04-145-11	Catalytic domain	1-398	NP_002740	N-terminal GST	Insect (sf21)
IKKα(CHUK) [inactive mutant]	BTN	BTN-IKKα(CHUK)[inactivemutant]	05-412-11-20N	Full-length	1-745	NP_001269.3	N-terminal DYKDDDDK	Insect (sf21)
JNK1(MAPK8) [inactive mutant]	GST	JNK1(MAPK8)[inactivemutant]	04-163-11	Catalytic domain	2-364	NP_620634.1	N-terminal GST	E. coli
MAP2K1 [inactive mutant]	GST	MAP2K1[inactivemutant]	07-109-11	Full-length	1-393	NP_002746.1	N-terminal GST	Insect (sf21)
MAP2K6 [inactive mutant]	GST	MAP2K6[inactivemutant]	07-146-11	Full-length	1-334	NP_002749.2	N-terminal GST	E. coli
MAP2K7 [inactive mutant]	GST	MAP2K7[inactivemutant]	07-110-11	Full-length	1-419	NP_660186.1	N-terminal GST	Insect (sf21)
p38α(MAPK14) [inactive mutant]	GST	p38α(MAPK14)[inactivemutant]	04-152-11	Truncated	9-352	NP_620581.1	N-terminal GST	E. coli

** Carna' s 'inactive' kinases are confirmed to be less active than our 'active' kinase products, and they can become fully catalytically active following phosphorylation by an upstream kinase.
In contrast, our 'inactive mutant' kinases are modified to be catalytically inactive, and as such are unable to phosphorylate substrates despite being phosphorylated by upstream kinases.

Protein Substrates

Protein Kinase	Tag	CarnaProductName	Catalog No.	Construct	Sequence	Accession Number	Tag Location	Expression System
Cofilin2	GST	Cofilin2	00-003	Full-length	1-166	NP_068733.1	N-terminal GST	E. coli
EIF2S1	GST	EIF2S1	00-006	Full-length	1-315	NP_004085.1	N-terminal GST	E. coli
EIF4EBP1	GST	EIF4EBP1	00-008	Full-length	1-118	NP_004086.1	N-terminal GST	E. coli
Smad1	GST	Smad1	00-000	Full-length	1-465	NP_005891.1	N-terminal GST	E. coli
Smad3	GST	Smad3	00-004	Full-length	1-425	NP_005893.1	N-terminal GST	E. coli

“QuickScout Screening Assist™ Kits” Kinase Protein Assay Kits

FP(IMAP™)

AKT1	AKT2	AKT3	AurA(AURKA)	AurB(AURKB)/INCENP
AurC(AURKC)	BRSK2	CaMK2α(CAMK2A)	CaMK4	CDK1(CDC2)/CycB1
CDK2/CycA2	CDK3/CycE1	CDK4/CycD3	CDK5/p25	CGK2(PRKG2)
CHK1(CHEK1)	CHK2(CHEK2)	CK1δ(CSNK1D)	CLK1	CRIK(CIT)
DAPK1	Erk1(MAPK3)	Erk2(MAPK1)	HGK(MAP4K4)	IKKα(CHUK)
IKKε(IKBKE)	IRAK1	IRAK4	MAPKAPK2	MAPKAPK3
MAPKAPK5	MELK	MNK1(MKNK1)	MNK2(MKNK2)	MRCKα(CDC42BPA)
MSK1(RPS6KA5)	MSK2(RPS6KA4)	MST4	NIM1K(MGC42105)	NuaK1
p38α(MAPK14)	p38β(MAPK11)	p38γ(MAPK12)	p38δ(MAPK13)	p70S6K(RPS6KB1)
p70S6Kβ(RPS6KB2)	PASK	PBK	PEK(EIF2AK3)	PGK(PRKG1)
PHKG1	PIM1	PIM2	PKACα(PRKACA)	PKCa(PRKCA)
PKCβ1(PRKCB1)	PKCδ(PRKCD)	PKCε(PRKCE)	PKCζ(PRKCZ)	PKCη(PRKCH)
PKCθ(PRKCQ)	PKCi(PRKCI)	PKR(EIF2AK2)	PLK1	PLK2
PLK3	PRKD1(PKD1)	PRKD2(PKD2)	PRKD3(PKD3)	PRKX
ROCK1	ROCK2	RSK1(RPS6KA1)	RSK2(RPS6KA3)	RSK3(RPS6KA2)
SGK	SGK3(SGKL)	SRPK1	SRPK2	TNIK
WNK1				

TR-FRET

ACTR2B (ACVR2B)	ACVR1B(ALK4)	ACVR2A	ACVRL1(ALK1)	ALK
ALK2(ACVR1)	AXL	BMPR1A	BMPR1B	BTK
CSK	EPHA2	FAK(PTK2)	FES	FGFR3
FLT3	HCK	IGF1R	JAK1	JAK2
JAK3	LCK	LYNa	MER(MERTK)	MET
PDGFRβ(PDGFRB)	SYK	TGFβR1(TGFBR1)	TIE2(TEK)	TRKA(NTRK1)
TYK2				

ELISA

BRAF	BRAF [V600E]	CDK8/CycC	CDK19(CDC2L6)/CycC	COT(MAP3K8)
DLK(MAP3K12)	Erk5(MAPK7)	GSK3β(GSK3B)	LZK(MAP3K13)	MAP2K1
MAP2K2	MAP2K3	MAP2K4	MAP2K5	MAP2K6
MAP2K7	MAP3K1	MAP3K2	MAP3K3	MAP3K4
MLK1(MAP3K9)	MLK2(MAP3K10)	MLK3(MAP3K11)	MOS	RAF1
TAK1-TAB1(MAP3K7)	TTK	WEE1		

ADP-Glo™

DGKα(DGKA)	DGKβ(DGKB)	DGKγ(DGKG)	DGKδ(DGKD)	DGKε(DGKE)
DGKη(DGKH)	DGKi(DGKI)	DGKθ(DGKQ)	DGKζ(DGKZ)	PIK3CA/PIK3R1
PIK3CA[E542K]/PIK3R1	PIK3CA[E545K]/PIK3R1	PIK3CA[H1047R]/PIK3R1	PIK3CA[P539R]/PIK3R1	PIK3CA[R88Q]/PIK3R1
PIK3CB/PIK3R1	PIK3CD/PIK3R1	PIKFYVE(PIP5K3)	PIP4K2A	PIP4K2B
PIP5K1A	PIP5K1B	PIP5K1C	PIP5KL1	

For details of services, please see "PRODUCTS AND SERVICES" on page 4.

Note: Please contact us for availability or further information. Information may be changed without notice.

“QuickScout Custom Profiling & Panel Profiling Series” Targets

Tyrosine Kinases	Custom		Panel
	Km bin	1mM	TK
ABL(ABL1)	●	●	●
ABL(ABL1) [E255K]	●	●	●
ABL(ABL1) [T315I]	●	●	●
ACK(TNK2)	●	●	●
ALK	●	●	●
ALK [C1156Y]	●	●	●
ALK [F1174L]	●	●	●
ALK [G1202R]	●	●	●
ALK [G1269A]	●	●	●
ALK [L1196M]	●	●	●
ALK [R1275Q]	●	●	●
ALK [T1151_L1152insT]	●	●	●
ARG(ABL2)	●	●	●
AXL	●	●	●
BLK	●	●	●
BMX	●	●	●
BRK(PTK6)	●	●	●
BTK	●	●	●
BTK [C481S]	●	●	●
CSK	●	●	●
DDR1	●	●	●
DDR2	●	●	●
EGFR	●	●	●
EGFR [C797S/L858R]	●	●	●
EGFR [d746-750]	●	●	●
EGFR [d746-750/C797S]	●	●	●
EGFR [d746-750/T790M]	●	●	●
EGFR [D770_N771insNPG]	●	●	●
EGFR [L858R]	●	●	●
EGFR [L861Q]	●	●	●
EGFR [T790M]	●	●	●
EGFR [T790M/C797S/L858R]	●	●	●
EGFR [T790M/L858R]	●	●	●
EML4-ALK	●	●	●
EPHA1	●	●	●
EPHA2	●	●	●
EPHA3	●	●	●
EPHA4	●	●	●
EPHA5	●	●	●
EPHA6	●	●	●
EPHA7	●	●	●
EPHA8	●	●	●
EPHB1	●	●	●
EPHB2	●	●	●
EPHB3	●	●	●
EPHB4	●	●	●
FAK(PTK2)	●	●	●
FER	●	●	●
FES	●	●	●
FGFR1	●	●	●
FGFR1 [V561M]	●	●	●
FGFR2	●	●	●
FGFR2 [V564I]	●	●	●
FGFR3	●	●	●
FGFR3 [K650E]	●	●	●
FGFR3 [K650M]	●	●	●
FGFR3 [V555L]	●	●	●
FGFR3 [V555M]	●	●	●
FGFR4	●	●	●
FGFR4 [N535K]	●	●	●
FGFR4 [V550E]	●	●	●
FGFR4 [V550L]	●	●	●
FGR	●	●	●
FLT1	●	●	●
FLT3	●	●	●
FLT4	●	●	●
FMS(CSF1R)	●	●	●
FRK	●	●	●
FYN [isoform a]	●	●	●
FYN [isoform b]	●	●	●
HCK	●	●	●
HER2(ERBB2)	●	●	●
HER4(ERBB4)	●	●	●
IGF1R	●	●	●
INSR	●	●	●
IRR(INSRR)	●	●	●
ITK	●	●	●
JAK1	●	●	●
JAK2	●	●	●
JAK3	●	●	●
KDR	●	●	●
KIT	●	●	●
KIT [D816E]	●	●	●
KIT [D816V]	●	●	●
KIT [D816Y]	●	●	●
KIT [T670I]	●	●	●
KIT [V560G]	●	●	●
KIT [V654A]	●	●	●
LCK	●	●	●
LTK	●	●	●
LYNa	●	●	●
LYNb	●	●	●
MER(MERTK)	●	●	●
MET	●	●	●
MET [D1228H]	●	●	●
MET [M1250T]	●	●	●
MET [Y1235D]	●	●	●
MUSK	●	●	●
NPM1-ALK	●	●	●
PDGFRa(PDGFR)	●	●	●
PDGFRa(PDGFR) [D842V]	●	●	●
PDGFRa(PDGFR) [T674I]	●	●	●
PDGFRa(PDGFR) [V561D]	●	●	●
PDGFRb(PDGFRB)	●	●	●
PYK2(PTK2B)	●	●	●
RET	●	●	●
RET [G691S]	●	●	●
RET [M918T]	●	●	●
RET [S891A]	●	●	●
RET [Y791F]	●	●	●
RON(MST1R)	●	●	●
ROS(ROS1)	●	●	●
SRC	●	●	●
SRM(SRMS)	●	●	●
SYK	●	●	●
TEC	●	●	●
TIE2(TEK)	●	●	●
TNK1	●	●	●
TRKA(NTRK1)	●	●	●
TRKB(NTRK2)	●	●	●

Tyrosine Kinases	Custom		Panel
	Km bin	1mM	TK
TRKC(NTRK3)	●	●	●
TXK	●	●	●
TYK2	●	●	●
TYRO3	●	●	●
YES(YES1)	●	●	●
YES(YES1) [T348I]	●	●	●
ZAP70	●	●	●

Serine/Threonine Kinases	Custom		Panel		
	Km bin	1mM	STK	MAPK	Cell
AKT1	●	●	●	●	●
AKT2	●	●	●	●	●
AKT3	●	●	●	●	●
AMPKa1/B1/v1(PRKAA1/B1/G1)	●	●	●	●	●
AMPKa2/B1/v1(PRKAA2/B1/G1)	●	●	●	●	●
AurA(AURKA)	●	●	●	●	●
AurA(AURKA)/TPX2	●	●	●	●	●
AurB(AURKB)/INCENP	●	●	●	●	●
AurC(AURKC)	●	●	●	●	●
BRAF	●	●	●	●	●
BRAF [V600E]	●	●	●	●	●
BRSK1	●	●	●	●	●
BRSK2	●	●	●	●	●
BUB1/BUB3	●	●	●	●	●
CaMK1a(CAMK1)*	●	●	●	●	●
CaMK1d(CAMK1D)	●	●	●	●	●
CaMK2a(CAMK2A)	●	●	●	●	●
CaMK2b(CAMK2B)	●	●	●	●	●
CaMK2γ(CAMK2G)	●	●	●	●	●
CaMK2δ(CAMK2D)	●	●	●	●	●
CaMK4	●	●	●	●	●
CDK7/ASK	●	●	●	●	●
CDK1(CDC2)/CycB1	●	●	●	●	●
CDK2/CycA2	●	●	●	●	●
CDK2/CycE1	●	●	●	●	●
CDK3/CycE1*	●	●	●	●	●
CDK4/CycD1*	●	●	●	●	●
CDK4/CycD3	●	●	●	●	●
CDK5/p25	●	●	●	●	●
CDK6/CycD3	●	●	●	●	●
CDK7/CyclH/MAT1	●	●	●	●	●
CDK9/CycT1	●	●	●	●	●
CGK2(PRKG2)	●	●	●	●	●
CHK1(CHEK1)	●	●	●	●	●
CHK2(CHEK2)	●	●	●	●	●
CK1a(CSNK1A1)	●	●	●	●	●
CK1γ1(CSNK1G1)	●	●	●	●	●
CK1γ2(CSNK1G2)	●	●	●	●	●
CK1γ3(CSNK1G3)	●	●	●	●	●
CK1δ(CSNK1D)	●	●	●	●	●
CK1ε(CSNK1E)	●	●	●	●	●
CK2a1/B(CSNK2A1/B)	●	●	●	●	●
CK2a2/B(CSNK2A2/B)	●	●	●	●	●
CLK1	●	●	●	●	●
CLK2	●	●	●	●	●
CLK3	●	●	●	●	●
COT(MAP3K8)	●	●	●	●	●
CRIK(CIT)	●	●	●	●	●
DAPK1	●	●	●	●	●
DCAMKL1	●	●	●	●	●
DCAMKL2	●	●	●	●	●
DLK(MAP3K12)	●	●	●	●	●
DYRK1A	●	●	●	●	●
DYRK1B	●	●	●	●	●
DYRK2	●	●	●	●	●
DYRK3	●	●	●	●	●
EEF2K	●	●	●	●	●
Erk1(MAPK3)	●	●	●	●	●
Erk2(MAPK1)	●	●	●	●	●
Erk5(MAPK7)*	●	●	●	●	●
GSK3a(GSK3A)	●	●	●	●	●
GSK3β(GSK3B)	●	●	●	●	●
Hsp90α(GSG2)	●	●	●	●	●
HGK(MAP4K4)	●	●	●	●	●
HIPK1	●	●	●	●	●
HIPK2	●	●	●	●	●
HIPK3	●	●	●	●	●
HIPK4	●	●	●	●	●
HPK1(MAP4K1)	●	●	●	●	●
IKKα(CHUK)	●	●	●	●	●
IKKβ(IKBKB)	●	●	●	●	●
IKKε(IKBKE)	●	●	●	●	●
IRAK1	●	●	●	●	●
IRAK4*	●	●	●	●	●
JNK1(MAPK8)	●	●	●	●	●
JNK2(MAPK9)	●	●	●	●	●
JNK3(MAPK10)	●	●	●	●	●
LATS1/MOBKL1A	●	●	●	●	●
LATS2/MOBKL1A	●	●	●	●	●
LOK(STK10)	●	●	●	●	●
MAP2K1	●	●	●	●	●
MAP2K2	●	●	●	●	●
MAP2K3	●	●	●	●	●
MAP2K4	●	●	●	●	●
MAP2K5	●	●	●	●	●
MAP2K6	●	●	●	●	●
MAP2K7	●	●	●	●	●
MAP3K1	●	●	●	●	●
MAP3K2	●	●	●	●	●
MAP3K3	●	●	●	●	●
MAP3K4	●	●	●	●	●
MAP3K5	●	●	●	●	●
MAP4K2	●	●	●	●	●
MAP4K5(KHS1)	●	●	●	●	●
MAPKAPK2	●	●	●	●	●
MAPKAPK3	●	●	●	●	●
MAPKAPK5	●	●	●	●	●
MARK1	●	●	●	●	●
MARK2	●	●	●	●	●
MARK3	●	●	●	●	●
MARK4	●	●	●	●	●
MELK	●	●	●	●	●
MINK(MINK1)	●	●	●	●	●

For details of services, please see "PRODUCTS AND SERVICES" on page 5~6. Note: Please contact us for availability or further information. Information may be changed without notice.

Serine/Threonine Kinases	Custom		Panel		
	Km bin	1mM	STK	MAPK	Cell
MLK1(MAP3K9)		●		●	
MLK2(MAP3K10)		●		●	
MLK3(MAP3K11)		●		●	
MNKK1(MKNK1)	●				
MNKK2(MKNK2)		●			
MOS		●		●	
MRCKα(CDC42BPB)	●				
MRCKβ(CDC42BPB)	●				
MSK1(RPS6KA5)		●			
MSK2(RPS6KA4)		●			
MSSK1(STK23)	●				
MST1(STK4)		●	●		
MST2(STK3)	●	●			
MST3(STK24)	●				
MST4	●				
NDR1(STK38)	●				
NDR2(STK38L)	●				
NEK1	●	●			●
NEK2	●	●	●		●
NEK4	●				
NEK6	●	●			●
NEK7	●	●			●
NEK9	●	●			●
NIM1K(MGC42105)	●				
NuaK1	●	●			
NuaK2	●	●			
p38α(MAPK14)	●	●	●		●
p38β(MAPK11)	●	●		●	
p38γ(MAPK12)	●	●		●	
p38δ(MAPK13)	●	●		●	
p70S6K(RPS6KB1)	●	●	●		●
p70S6Kβ(RPS6KB2)	●				
PAK1	●	●			
PAK2	●		●		
PAK4	●				
PAK5(PAK7)	●	●			
PAK6	●				
PASK	●	●			
PBK	●	●	●		
PDHK2(PDK2)	●				
PDHK4(PDK4)	●				
PDK1(PDPK1)	●	●	●		
PEK(EIF2AK3)	●				
PGK(PRKG1)	●				
PHKG1	●				
PHKG2	●				
PIM1	●	●	●		●
PIM2	●				
PIM3	●		●		●
PKACα(PRKACA)	●	●	●		
PKACβ(PRKACB)	●	●			
PKACγ(PRKACG)	●				
PKCα(PRKCA)	●	●	●		
PKCβ1(PRKCB1)	●				
PKCβ2(PRKCB2)	●				
PKCγ(PRKCG)	●	●			
PKCδ(PRKCD)	●				
PKCε(PRKCE)	●	●			
PKCζ(PRKCZ)	●				
PKCη(PRKCH)	●				
PKCθ(PRKCO)	●				
PKCι(PRKCI)	●				
PKN1	●				
PKR(EIF2AK2)	●				
PLK1	●	●			●
PLK2	●				●
PLK3	●	●			
PRKD1(PKD1)	●		●		
PRKD2(PKD2)	●	●			
PRKD3(PKD3)	●				
PRKX	●				
QIK(SNF1LK2)	●	●			
RAF1	●			●	
ROCK1	●	●	●		
ROCK2	●	●			
RSK1(RPS6KA1)	●	●			●
RSK2(RPS6KA3)	●	●			
RSK3(RPS6KA2)	●	●			
RSK4(RPS6KA6)	●	●			
SGK	●	●	●		
SGK2	●				
SGK3(SGKL)	●				
SIK(SNF1LK)	●	●			
SIK3(OSK)	●	●			
skMLCK(MYLK2)*	●	●			
SLK	●				
SRPK1	●				
SRPK2	●				
TAK1-TAB1(MAP3K7)	●	●			
TAOK2	●				
TBK1	●	●			
TNKK	●	●			
TSSK1B(TSSK1)	●	●	●		
TSSK2	●				
TSSK3	●				
WNK1	●				
WNK2	●				
WNK3	●				

“QuickScout Custom Profiling Series” Preincubation Targets

Tyrosine Kinases
ABL(ABL1)
ABL(ABL1) [E255K]
ABL(ABL1) [T315I]
ALK
ALK [C1156Y]
ALK [G1202R]
ALK [G1269A]
ALK [F1174L]
ALK [L1196M]
ALK [R1275Q]
ARG(ABL2)
BMX
EGFR
EGFR [d746-750/C797S]
EGFR [T790M/L858R]
EPHA1
EPHA2
EPHA3
EPHA4
EPHA5
EPHA6
EPHA7
EPHA8
EPHB1
EPHB2
EPHB3
EPHB4
FER
FES
FGFR1
FGFR2 [V564I]
FGFR4
FGFR4 [V550L]
FLT1
FLT4
FRK
HER4(ERBB4)
INSR
JAK2
KDR
LCK
LYNa
LYNb
MER(MERTK)
MET
MET [D1228H]
MET [M1250T]
MET [Y1235D]
NPM1-ALK
PDGFRα(PDGFRα)
PDGFRα(PDGFRα) [D842V]
PDGFRβ(PDGFRβ)
PYK2(PTK2B)

Tyrosine Kinases
RET
RET [G691S]
RET [M918T]
RET [S891A]
RET [Y791F]
RON(MST1R)
ROS(ROS1)
SRC
TEC
TIE2(TEK)
TRKB(NTRK2)
TRKC(NTRK3)
TYRO3
YES(YES1)

Serine/Threonine Kinases
AKT1
AKT2
AKT3
AMPA1/β1/γ1(PRKAA1/B1/G1)
AMPA2/β1/γ1(PRKAA2/B1/G1)
BRK1
BRK2
BUB1/BUB3
CaMK2α(CAMK2A)
CaMK2β(CAMK2B)
CaMK2γ(CAMK2G)
CaMK2δ(CAMK2D)
CaMK4
CDK1(CDC2)/CycB1
CDK2/CycA2
CDK2/CycE1
CDK5/p25
CGK2(PRKG2)
CHK1(CHEK1)
CHK2(CHEK2)
CK1γ1(CSNK1G1)
CK1γ2(CSNK1G2)
CK1γ3(CSNK1G3)
CK1δ(CSNK1D)
CK2α1/β(CSNK2A1/B)
CK2α2/β(CSNK2A2/B)
CLK1
CLK2
CLK3
DYRK1A
DYRK1B
DYRK2
DYRK3
Erk1(MAPK3)
Erk2(MAPK1)

Serine/Threonine Kinases
Haspin(GSG2)
HIPK1
HIPK2
HIPK3
IKKβ(IKBKB)
JNK1(MAPK8)
JNK2(MAPK9)
JNK3(MAPK10)
MAP4K5(KHS1)
MAPKAPK3
MAPKAPK5
MARK1
MARK3
MARK4
MNK1(MKNK1)
MNK2(MKNK2)
MSK1(RPS6KA5)
NEK2
NEK4
NIM1K(MGC42105)
Nuak1
Nuak2
p38α(MAPK14)
p38β(MAPK11)
p38γ(MAPK12)
p38δ(MAPK13)
p70S6K(RPS6KB1)
p70S6Kβ(RPS6KB2)
PAK1
PAK2
PAK5(PAK7)
PIM1
PKCε(PRKCε)
PRKD1(PKD1)
PRKD2(PKD2)
QIK(SNF1LK2)
SGK
SGK3(SGKL)
skMLCK(MYLK2)
TBK1
TSSK1B(TSSK1)

Tyrosine Kinases 67

Serine/Threonine Kinases 76

Total 143 targets

For details of services, please see "PRODUCTS AND SERVICES" on page 5.

Note: Please contact us for availability or further information.

NanoBRET™ TE Intracellular Kinase Cell-Based Assay Service Targets

Targets in blue letters have been successfully validated at Carna.
Please contact us about the turnaround time for the targets in grey letters.

Kinase Name					
AAK1	CDK7	FAK(PTK2)	LIMK2	NLK	RSK2(RPS6KA3) [L608F]
ABL1(ABL)	CDK7/Cyc H	FER	LRRK2	NRK	RSK3(RPS6KA2)
ABL1(ABL) [E255K]	CDK8/Cyc C	FES	LRRK2 [G2019S]	NUAK1(ARK5)	RSK4(RPS6KA6)
ABL1(ABL) [F317I]	CDK9/Cyc K	FGFR1	LRRK2 [I2020T]	NUAK2	SBK3
ABL1(ABL) [F317L]	CDK9/Cyc T1	FGFR2	LRRK2 [R1441C]	PAK4	SGK1(SGK)
ABL1(ABL) [H396P]	CDK10/Cyc L2	FGFR2 [K659M]	LTK	PAK6	SGK2
ABL1(ABL) [M351T]	CDK11A/Cyc K	FGFR2 [L617F]	LYN	PAK7(PAK5)	SIK1(SIK)
ABL1(ABL) [Q252H]	CDK11A/Cyc L2	FGFR2 [L617V]	MAP2K5(MEK5)	PDGFRA [V561D]	SIK2(QIK, SNF1LK2)
ABL1(ABL) [T315I]	CDK12/Cyc K	FGFR2 [M537I]	MAP2K6	PHKG1	SIK3(QSK)
ABL1(ABL) [Y253F]	CDK13/Cyc K	FGFR2 [N549H]	MAP3K2	PHKG2	SLK
ABL2(ARG)	CDK14/Cyc Y	FGFR2 [N549K]	MAP3K3	PI4KA	SNRK
ACVR1(ALK2)	CDK15/Cyc Y	FGFR2 [V564F]	MAP3K4	PI4KB	SRC
ACVR1(ALK2) [G328V]	CDK16(PCTAIRE1)/CycY	FGFR2 [V564I]	MAP3K9(MLK1)	PIK3C3	SRMS(SRM)
ACVR1(ALK2) [G356D]	CDK17/Cyc Y	FGFR3	MAP3K10(MLK2)	PIK3CA/PIK3R1	STK3(MST2)
ACVR1(ALK2) [Q207D]	CDK18/Cyc Y	FGFR3 [G380R]	MAP3K11(MLK3)	PIK3CA [C420R]/PIK3R1	STK4(MST1)
ACVR1(ALK2) [R206H]	CDK19/Cyc C	FGFR3 [G697C]	MAP3K12(DLK)	PIK3CA [E542K]/PIK3R1	STK10(LOK)
ACVRL1(ALK1)	CDK20/Cyc H	FGFR3 [V555M]	MAP3K13(LZK)	PIK3CA [E545A]/PIK3R1	STK11
ADK	CDKL1	FGFR4	MAP3K19	PIK3CA [E545K]/PIK3R1	STK16
AKT1	CDKL2	FGR	MAP3K21(MLK4)	PIK3CA [H1047L]/PIK3R1	STK17A(DRAK1)
AKT1 [E17K]	CDKL3	FLT1	MAP4K1(HPK1)	PIK3CA [H1047R]/PIK3R1	STK17B
AKT2	CDKL5	FLT3	MAP4K2	PIK3CA [H1047Y]/PIK3R1	STK24(MST3)
AKT2 [E17K]	CHEK1(CHK1)	FLT3 [D835H]	MAP4K3	PIK3CA [I800L]/PIK3R1	STK26(MST4)
AKT3 [E17K]	CHEK2(CHK2)	FLT3 [D835V]	MAP4K5	PIK3CA [M1043I]/PIK3R1	STK32A
AKT3 [G171R]	CLK1	FLT3 [D835Y]	MAPK4	PIK3CA [Q546K]/PIK3R1	STK32B
ALK	CLK2	FLT3 [K663Q]	MAPK6	PIK3CB/PIK3R1	STK33
ALK4(ACVR1B)	CLK4	FLT3 [N841I]	MAPK11(p38β)	PIK3CD/PIK3R1	STK35
AURKA	COQ8B	FLT3 [R834Q]	MAPK14(p38α)	PIK3CG	STK36
AURKB	CRAF(RAF1)	FRK	MAPK14(p38α) [T106M]	PIKFYVE(PIP5K3)	STK38(NDR1)
AURKC	CSF1R(FMS)	FYN	MARK2	PIM3	STK38L(NDR2)
AXL	CSK	FYN [Y531F]	MARK3	PIP4K2C	TBK1
BLK	CSNK1A1L	GAK	MARK4	PIP5K1B	TEC
BMP2K	CSNK1D(CK1δ)	GSK3A(GSK3α)	MAST3	PKACα(PRKACA)	TEK(TIE2)
BMPRI1A	CSNK1E(CK1ε)	GSK3B(GSK3β)	MAST4	PKACβ(PRKACB)	TEK(TIE2) [A1124V]
BMX	CSNK1G2(CK1γ2)	HCK	MELK	PKCα(PRKCA)	TEK(TIE2) [P883A]
BRAF	CSNK2A1(CK2α1)	HIPK2	MELK [T460M]	PKCβ1(PRKCB1)	TEK(TIE2) [R849W]
BRAF [V600E]	CSNK2A2(CK2α2)	HIPK3	MER(MERTK)	PKCγ(PRKCG)	TEK(TIE2) [Y1108F]
BRK(PTK6)	DAPK2	HIPK4	MER(MERTK) [A708S]	PKCδ(PRKCD)	TEK(TIE2) [Y897C]
BRSK1	DCLK1(DCAMKL1) [isoform1]	ICK	MET	PKCε(PRKCE)	TEK(TIE2) [Y897S]
BRSK2	DCLK1(DCAMKL1) [isoform2]	IGF1R	MET [D1228H]	PKCη(PRKCH)	TESK1
BTk	DCLK3	IKBKE(IKKe)	MET [D1228N]	PKCθ(PRKCQ)	TGFBR1(ALK5)
BTk [C481S]	DDR1	INSR	MET [F1200I]	PKMYT1(MYT1)	TGFBR2
BTk [E41K]	DDR2	IRAK1	MET [M1250T]	PKN1	TIE1
BTk [L528W]	DDR2 [N456S]	IRAK3	MET [P991S]	PLK1	TLK1
BTk [P190K]	DYRK1A	IRAK4	MET [T1173I]	PLK2	TLK2
BTk [T474I]	DYRK1B	ITK	MET [T992I]	PLK3	TNK1
CAMK1(CaMK1α)	DYRK2	JAK1	MET [V1092I]	PLK4	TNK2(ACK)
CAMK1D(CaMK1δ)	EGFR	JAK1_JH2 Domain	MET [Y1230A]	PNCK(CaMK1β)	TNNI3K(HH498)
CAMK1G(CaMK1γ)	EGFR [d746-750]	JAK2	MET [Y1230C]	PRKAA1(AMPKα1)	TRKA(NTRK1)
CAMK2A(CaMK2α)	EGFR [L858R]	JAK2 [V617F]	MET [Y1230D]	PRKAA2(AMPKα2)	TRKA(NTRK1) [G667C]
CAMK2D(CaMK2δ)	EGFR [T790M/L858R]	JAK2_JH1 Domain	MET [Y1230H]	PRKG2(CGK2)	TRKB(NTRK2)
CAMK2G(CaMK2γ)	EIF2AK4(GCN2)_Domain2	JAK3	MET [Y1235D]	PRKX	TRKC(NTRK3)
CAMKK1(124-411)	EPHA1	JNK1(MAPK8)	MKNK2(MNK2)	PYK2 (PTK2B)	TSSK1B
CAMKK2(165-445)	EPHA2	JNK2(MAPK9)	MLTK(ZAK)	RET	TTK
CDK1/Cyc A2(CDC2/CycA2)	EPHA3	JNK3(MAPK10)	MOK	RET [M918T]	TXK
CDK1/Cyc B1(CDC2/CycB1)	EPHA4	KDR(VEGFR2)	MUSK	RET [V804L]	TYK2
CDK1/Cyc E1(CDC2/CycE1)	EPHA5	KIT	MYLK2(skMLCK)	RET [V804M]	TYK2_JH1 Domain
CDK2/Cyc A1	EPHA6	KIT [A829P]	MYLK3	RIOK2	TYK2_JH2 Domain
CDK2/Cyc A2	EPHA7	KIT [D816H]	MYLK4	RIPK1	TYRO3
CDK2/Cyc E1	EPHA8	KIT [D816V]	NEK1	RIPK2	ULK1
CDK3/Cyc E1	EPHB1	KIT [L576P]	NEK2	RIPK3	ULK2
CDK4/Cyc D1	EPHB2	KIT [V559D]	NEK3	ROCK1	ULK3
CDK4/Cyc D3	EPHB3	KIT [V559D,T670I]	NEK4	ROCK2	WEE1
CDK5	EPHB4	KIT [V559D,V654A]	NEK5	RON	WEE2
CDK5/CDK5R1	ERK1(MAPK3)	LATS1	NEK6	RPS6KA4(MSK2)	YES1(YES)
CDK5/CDK5R2	ERK2(MAPK1)	LATS2	NEK9	RSK1(RPS6KA1)	
CDK6/Cyc D1	ERN1	LCK	NEK11	RSK2(RPS6KA3)	
CDK6/Cyc D3	ERN2	LIMK1	NIM1K(MGC42105)	RSK2(RPS6KA3) [I416V]	

For details of services, please see "PRODUCTS AND SERVICES" on page 7.
Note: Please contact us for availability or further information.

NanoBRET™ TE Intracellular Kinase Cell-Based Assay Service Targets 【Panel Services】

～CDK Panel Service～

Kinase Name			
CDK1(CDC2)／CycB1	CDK5／CDK5R1	CDK9／CycT1	CDK15／CycY
CDK2／CycA1	CDK6／CycD1	CDK10／CycL2	CDK16／CycY
CDK2／CycE1	CDK6／CycD3	CDK11A／CycL2	CDK17／CycY
CDK3／CycE1	CDK7／CycH	CDK12／CycK	CDK18／CycY
CDK4／CycD1	CDK8／CycC	CDK13／CycK	CDK19／CycC
CDK4／CycD3	CDK9／CycK	CDK14／CycY	CDK20／CycH

～ Kinome-Wide Profiling Service (K192) ～

Kinase Name			
AAK1	DCLK3	LTK	PHKG2
ACK (TNK2)	DLK (MAP3K12)	MAP3K2	PKACα (PRKACA)
AKT2	DRAK2 (STK17B)	MAP3K3	PKACβ (PRKACB)
AMPKα1 (PRKAA1)	DYRK1A	MAP3K4	PKCε (PRKCE)
AMPKα2 (PRKAA2)	DYRK1B	MAP3K19	PLK2
ARG (ABL2)	EPHA1	MAP4K2	PLK3
AurA (Aurora A)	EPHA4	MAP4K3	PLK4
AurB (Aurora B)	EPHA6	MAP4K5	PRKX
AurC (Aurora C)	EPHA7	MAPK4	PYK2 (PTK2B)
AXL	EPHB1	MAPK6	QIK (SIK2)
BMP2K	EPHB4	MARK2	RET
BMX	ERN1	MARK4	RIOK2
BRAF[V600E]	ERN2	MAST3	RIPK1
BRK (PTK6)	FAK (PTK2)	MAST4	RIPK2
BRSK1	FER	MELK	RON
BRSK2	FES	MER (MERTK)	RSK1
BTk	FGFR1	MET	RSK2
CaMK1α (CAMK1)	FGFR2	MLK1 (MAP3K9)	RSK3
CaMK2α (CAMK2A)	FGFR3	MLK2 (MAP3K10)	RSK4
CaMK2δ (CAMK2D)	FGFR4	MLK3 (MAP3K11)	SBK3
CDC2 (CDK1)／CycB1	FLT3	MLK4 (MAP3K21)	SGK (SGK1)
CDK2／CycE1	FYN [isoform a] (FYN)	MNK2 (MKNK2)	SIK (SIK1)
CDK3／CycE1	GAK	MPSK1 (STK16)	SIK3 (QSK)
CDK4／CycD3	HH498 (TNNI3K)	MSK2 (RPS6KA4)	skMLCK (MYLK2)
CDK5／p25 (CDK5R1)	HIPK2	MST1 (STK4)	SLK
CDK6／CycD1	HIPK3	MST2 (STK3)	SNRK
CDK7	HIPK4	MUSK	SRM (SRMS)
CDK9／CycK	HPK1 (MAP4K1)	MYLK3	STK32B
CDK10／CycL2	ICK	MYLK4	STK33
CDK14／CycY	IGF1R	MYT1 (PKMYT1)	STK35
CDK15／CycY	IKKε (IKBKE)	NDR1 (STK38)	STK36
CDK17／CycY	INSR	NDR2 (STK38L)	TBK1
CDK18／CycY	IRAK3	NEK1	TEC
CDK20／CycH	IRAK4	NEK2	TIE1
CDKL1	ITK	NEK3	TIE2 (TEK)
CDKL2	JAK2[V617F]	NEK4	TLK1
CDKL3	JAK3	NEK5	TLK2
CDKL5	JNK1	NEK9	TNK1
CHK2 (CHEK2)	JNK2	NEK11	TRKA (NTRK1)
CK1γ2 (CSNK1G2)	JNK3	NIM1K (MGC42105)	TRKB (NTRK2)
CK1δ (CSNK1D)	LATS1	NLK	TXK
CK2α1 (CSNK2A1)	LATS2	NUAK1	TYRO3
CK2α2 (CSNK2A2)	LCK	p38α (MAPK14)	ULK1
CLK1	LIMK1	p38β (MAPK11)	ULK2
CLK2	LIMK2	PAK4	ULK3
CLK4	LKB1 (STK11)	PAK6	WEE1
CSNK1A1L	LOK (STK10)	PCTAIRE1 (CDK16)／CycY	WEE2
DAPK2	LRRK2	PHKG1	ZAK (MLTK)

For details of services, please see "PRODUCTS AND SERVICES" on page 8.

Note: Please contact us for availability or further information.

Detection of Protein-Protein Interactions ~ProbeX™~ Stable Cell Lines List

GPCR

 A video of real time monitoring is available for the product with this mark on our website.

Cat.#	Target GPCR and Description	Receptor Family	Coupling	Accession Number	β-Arrestin Isoform	Cell Line	Species
PXC-001-1	ADCYAP1R1  Adenylate cyclase activating polypeptide 1 (pituitary) receptor type I	VIP and PACAP	Gs	NM_001118	β- Arrestin-1	HEK293	Human
PXC-002-2	ADRA2A  Alpha-2-adrenergic receptor	Adrenoceptor	Gi	NM_000681	β- Arrestin-2	HEK293	Human
PXC-003-1	ADRB2  Beta-2-adrenergic receptor	Adrenoceptor	Gs	NM_000024	β- Arrestin-1	HEK293	Human
PXC-003-2	ADRB2  Beta-2-adrenergic receptor	Adrenoceptor	Gs	NM_000024	β- Arrestin-2	HEK293	Human
PXC-004-1	AGTR1 Angiotensin II receptor, type 1	Angiotensin	Gq	NM_000685	β- Arrestin-1	HEK293	Human
PXC-004-2	AGTR1 Angiotensin II receptor, type 1	Angiotensin	Gq	NM_000685	β- Arrestin-2	HEK293	Human
PXC-005-1	AGTRL1 Apelin receptor	Apelin	Gi	NM_005161	β- Arrestin-1	HEK293	Human
PXC-005-2	AGTRL1 Apelin receptor	Apelin	Gi	NM_005161	β- Arrestin-2	HEK293	Human
PXC-006-1	CCKBR  Cholecystokinin B receptor	Cholecystokinin	Gq	NM_176875	β- Arrestin-1	HEK293	Human
PXC-006-2	CCKBR Cholecystokinin B receptor	Cholecystokinin	Gq	NM_176875	β- Arrestin-2	HEK293	Human
PXC-007-1	CCR7 CC chemokine receptor 7 (human)	Chemokine	Gi	NM_001838	β- Arrestin-1	HEK293	Human
PXC-007-2	CCR7 CC chemokine receptor 7 (human)	Chemokine	Gi	NM_001838	β- Arrestin-2	HEK293	Human
PXC-008-1	EDG1 Endothelial differentiation protein/sphingosine-1-phosphate receptor 1	Lysophospholipid	Gi	NM_001400	β- Arrestin-1	HEK293	Human
PXC-008-2	EDG1 Endothelial differentiation protein/sphingosine-1-phosphate receptor 1	Lysophospholipid	Gi	NM_001400	β- Arrestin-2	HEK293	Human
PXC-009-2	ENDRA Endothelin receptor type A	Endothelin	Gq	NM_001957	β- Arrestin-2	HEK293	Human
PXC-010-1	ENDRB Endothelin receptor type B	Endothelin	Gs	NM_000115	β- Arrestin-1	HEK293	Human
PXC-010-2	ENDRB Endothelin receptor type B	Endothelin	Gs	NM_000115	β- Arrestin-2	HEK293	Human
PXC-011-1	EP4  Prostaglandin E2 receptor 4	Prostanoid	Gs	NM_000958	β- Arrestin-1	HEK293	Human
PXC-011-2	EP4  Prostaglandin E2 receptor 4	Prostanoid	Gs	NM_000958	β- Arrestin-2	HEK293	Human
PXC-012-2	NPY1R Neuropeptide Y receptor Y1	Neuropeptide Y	Gi	NM_000909	β- Arrestin-2	HEK293	Human
PXC-013-1	OPRM1  Opioid receptor, mu 1	Opioid	Gi	NM_000914	β- Arrestin-1	HEK293	Human
PXC-013-2	OPRM1  Opioid receptor, mu 1	Opioid	Gi	NM_000914	β- Arrestin-2	HEK293	Human
PXC-014-1	SSTR2 Somatostatin receptor 2	Somatostatin	Gi	NM_001050	β- Arrestin-1	HEK293	Human
PXC-014-2	SSTR2 Somatostatin receptor 2	Somatostatin	Gi	NM_001050	β- Arrestin-2	HEK293	Human
PXC-015-1	EDG3 Endothelial differentiation protein/sphingosine-1-phosphate receptor 3	Lysophospholipid	Gi	NM_005226.2	β-Arrestin-1	HEK293	Human

Protein-Protein Interaction

Cat.#	Target Description	Accession Number (α)	Accession Number (β)	Cell Line	Species
PXC-PPI-1	Hif-1 complex Hypoxia inducible factor 1 alpha, Hydrocarbon receptor nuclear translocator (ARNT)	NM_001530	NM_001668	HEK293	Human
PXC-PPI-2	IκBα/b-TRCP Complex NFKB inhibitor alpha (NFKBIA), beta-transducin repeat containing E3 ubiquitin protein ligase (BTRC), transcript	NM_020529	NM_001256856	HEK293	Human
PXC-PPI-3	ARIA/PTEN Complex endothelial cell-specific molecule-2 (ECSM2), protein tyrosine phosphatase (TEP1)	DQ462572	U96180	HEK293	Human

Reagent

Cat.#	Name of Product
PXR-SG001	Split Glow Cell Assay Reagent

Split Glow will be included with each purchase of a stable cell line.

For details of services, please see "PRODUCTS AND SERVICES" on page 9.

Note: Please contact us for availability or further information.

Memo

[illegible]

Memo

[illegible]

Memo

[illegible]

Memo

[illegible]

Carna Biosciences, Inc.

BMA 3F, 1-5-5 Minatojima-Minamimachi,
Chuo-ku, Kobe 650-0047 JAPAN

TEL: +81 78 302 7091 / FAX: +81 78 302 7086

E-mail: info@carnabio.com