

Biotinylated Kinases

65
Available!

See reverse side for available kinases!

Use our products to study kinase-molecule interactions!

The immobilization of target proteins onto sensor surfaces, while maintaining both structure and activity, has been problematic and rate-limiting to the acquisition of meaningful SPR and/or BLI data in small molecule drug discovery. Carna Biosciences has developed proprietary technology to produce biotinylated kinase products to help overcome these challenges. A few features of our new and expanding product line are shown below.

Biotinylated Kinases

The Carna Advantage...

- Singly Biotinylated at N-Terminus
- Preserved kinase activity and structure
- Produced in insect cells using baculovirus expression system
- Highly purified by affinity column chromatography



Example of Application of *Biotinylated* Kinase (FLT3) in SPR

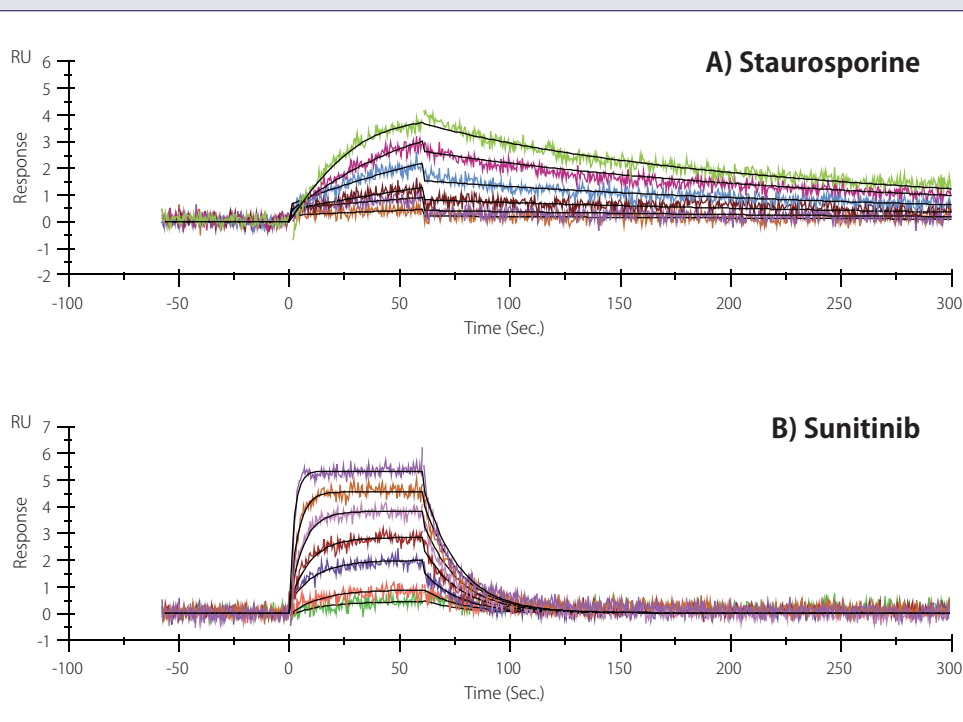


Fig. Representative SPR sensorgrams of the binding between biotinylated-FLT3 and kinase inhibitors. A) Staurosporine, and B) Sunitinib.

Production Steps

Cloning and Cell Culture

Purification

- *Standard Production*
- *ATP-Treated Kinases*
- *Phosphatase Treated Kinases*

will be offered on a fee-for-service basis in the future. Please contact us for more details.

Quality Control

Immediate Availability

Vialled as 10ug and 100ug aliquots

Purchase for any binding assay that utilizes biotin-avidin interactions!

For more details, please visit our website.

Carna Biosciences, Inc.

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

TEL: 078-302-7091 / FAX: 078-302-7086

E-mail: info@carnabio.com

Biotinylated Kinases from Carna Biosciences, Inc.

Cat No.	Name of Product	Length	Amino Acid	Accession Number	Tag	M.W.	Expression System
08-401-20N	BTN - ABL(ABL1)	Full-length	2-1130(end)	NP_005148.2	N-terminal DYKDDDDK	126 kDa	Insect (<i>Sf21</i>)
09-432-20N	BTN - ACTR2B(ACVR2B)	Cytoplasmic domain	161-512(end)	NP_001097.2	N-terminal DYKDDDDK	43 kDa	Insect (<i>Sf21</i>)
01-401-20N	BTN - AKT1	Catalytic domain	104-480(end)	NP_005154.1	N-terminal DYKDDDDK	47 kDa	Insect (<i>Sf21</i>)
01-402-20N	BTN - AKT2	Catalytic domain	120-481(end)	NP_001617.1	N-terminal DYKDDDDK	45 kDa	Insect (<i>Sf21</i>)
08-405-20N	BTN - ALK	Cytoplasmic domain	1058-1620(end)	BAG10812.1	N-terminal DYKDDDDK	67 kDa	Insect (<i>Sf21</i>)
08-429-20N	BTN - ALK[L1196M]	Cytoplasmic domain	1058-1620(end)	BAG10812.1	N-terminal DYKDDDDK	67 kDa	Insect (<i>Sf21</i>)
05-401-20N	BTN - AurA(AURKA)	Full-length	1-403(end)	NP_940835.1	N-terminal DYKDDDDK	50 kDa	Insect (<i>Sf21</i>)
05-402-21N	BTN - AurB(AURKB)/INCENP	Full-length	1-344	Q96GD4	N-terminal DYKDDDDK	43 kDa	Insect (<i>Sf21</i>)
08-407-20N	BTN - AXL	Cytoplasmic domain	464-885(end)	NP_001690.2	N-terminal DYKDDDDK	51 kDa	Insect (<i>Sf21</i>)
08-479-20N	BTN - BMX	Full-length	1-675(end)	NP_001712.1	N-terminal DYKDDDDK	82 kDa	Insect (<i>Sf21</i>)
08-480-20N	BTN - BTK	Full-length	2-659(end)	NP_000052.1	N-terminal DYKDDDDK	80 kDa	Insect (<i>Sf21</i>)
02-417-20N	BTN - CHK1(CHEK1)	Full-length	1-476(end)	NP_001265.1	N-terminal DYKDDDDK	58 kDa	Insect (<i>Sf21</i>)
08-414-20N	BTN - DDR2	Cytoplasmic domain	422-855(end)	NP_006173.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>Sf21</i>)
09-411-20N	BTN - DLK(MAP3K12)	Catalytic domain	1-520	NP_006292.3	N-terminal DYKDDDDK	63 kDa	Insect (<i>Sf21</i>)
08-415-20N	BTN - EGFR	Cytoplasmic domain	669-1210(end)	NP_005219.2	N-terminal DYKDDDDK	66 kDa	Insect (<i>Sf21</i>)
08-410-20N	BTN - EGFR[T790M/L858R]	Cytoplasmic domain	669-1210(end)	NP_005219.2	N-terminal DYKDDDDK	66 kDa	Insect (<i>Sf21</i>)
08-437-20N	BTN - FAK(PTK2)	Truncated protein	376-1052(end)	NP_722560.1	N-terminal DYKDDDDK	80 kDa	Insect (<i>Sf21</i>)
08-435-20N	BTN - FGFR1	Cytoplasmic domain	396-820(end)	NP_056934.2	N-terminal DYKDDDDK	52 kDa	Insect (<i>Sf21</i>)
08-433-20N	BTN - FGFR3	Cytoplasmic domain	436-806(end)	NP_000133.1	N-terminal DYKDDDDK	52 kDa	Insect (<i>Sf21</i>)
08-436-20N	BTN - FGFR4	Cytoplasmic domain	460-802(end)	NP_002002.3	N-terminal DYKDDDDK	42 kDa	Insect (<i>Sf21</i>)
08-454-20N	BTN - FLT3	Cytoplasmic domain	564-993(end)	NP_004110.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>Sf21</i>)
08-455-20N	BTN - FMS(CSF1R)	Cytoplasmic domain	538-972(end)	NP_005202.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>Sf21</i>)
08-468-20N	BTN - FYN[isoform a]	Full-length	1-537(end)	NP_002028.1	N-terminal DYKDDDDK	65 kDa	Insect (<i>Sf21</i>)
08-441-20N	BTN - IGF1R	Cytoplasmic domain	959-1367(end)	NP_000866.1	N-terminal DYKDDDDK	50 kDa	Insect (<i>Sf21</i>)
05-484-20N	BTN - IKK β (IKKB)	Truncated protein	1-662	NP_001547.1	N-terminal DYKDDDDK	80 kDa	Insect (<i>Sf21</i>)
05-412-20N	BTN - IKK α (CHUK)	Full-length	1-745(end)	NP_001269.3	N-terminal DYKDDDDK	88 kDa	Insect (<i>Sf21</i>)
05-412-11-20N	BTN - IKK α (CHUK)[inactive mutant]	Full-length	1-745(end)	NP_001269.3	N-terminal DYKDDDDK	88 kDa	Insect (<i>Sf21</i>)
09-445-20N	BTN - IRAK4	Full-length	1-460(end)	NP_057207.2	N-terminal DYKDDDDK	55 kDa	Insect (<i>Sf21</i>)
08-481-20N	BTN - ITK	Full-length	2-620(end)	NP_005537.3	N-terminal DYKDDDDK	75 kDa	Insect (<i>Sf21</i>)
08-444-20N	BTN - JAK1	Catalytic domain	850-1154(end)	NP_002218.2	N-terminal DYKDDDDK	39 kDa	Insect (<i>Sf21</i>)
08-445-20N	BTN - JAK2	Catalytic domain	826-1132(end)	NP_004963.1	N-terminal DYKDDDDK	39 kDa	Insect (<i>Sf21</i>)
08-491-20N	BTN - KDR(VEGFR2)	Cytoplasmic domain	790-1356(end)	NP_002244.1	N-terminal DYKDDDDK	67 kDa	Insect (<i>Sf21</i>)
08-456-21N	BTN - KIT	Cytoplasmic domain	544-976(end)	NP_000213	N-terminal DYKDDDDK	53 kDa	Insect (<i>Sf21</i>)
08-470-20N	BTN - LCK	Full-length	1-509(end)	NP_005347.2	N-terminal DYKDDDDK	62 kDa	Insect (<i>Sf21</i>)
08-471-20N	BTN - LYNa	Full-length	1-512(end)	NP_002341.1	N-terminal DYKDDDDK	63 kDa	Insect (<i>Sf21</i>)
07-441-10-20N	BTN - MAP2K1[inactive]	Full-length	1-393(end)	NP_002746.1	N-terminal DYKDDDDK	47 kDa	Insect (<i>Sf21</i>)
07-407-20N	BTN - MAP3K5	Catalytic domain	654-971	NP_005914.1	N-terminal DYKDDDDK	39 kDa	Insect (<i>Sf21</i>)
08-451-20N	BTN - MET	Cytoplasmic domain	956-1390(end)	NP_000236.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>Sf21</i>)
02-445-20N	BTN - MNK1(MKNK1)	Full-length	1-424(end)	BAA19885.1	N-terminal DYKDDDDK	51 kDa	Insect (<i>Sf21</i>)
05-433-20N	BTN - NFKB1	Truncated protein	1-346, 733-979(end)	NP_149107.4	N-terminal DYKDDDDK	70 kDa	Insect (<i>Sf21</i>)
04-452-20N	BTN - p38 α (MAPK14)	Truncated protein	9-352, 1-334(end)	NP_620581.1	N-terminal DYKDDDDK	43 kDa	Insect (<i>Sf21</i>)
04-452-10-20N	BTN - p38 α (MAPK14)[inactive]	Truncated protein	9-352	NP_620581.1	N-terminal DYKDDDDK	43 kDa	Insect (<i>Sf21</i>)
07-426-20N	BTN - PAK4	Full-length	1-591(end)	NP_005875.1	N-terminal DYKDDDDK	68 kDa	Insect (<i>Sf21</i>)
11-401-20N	BTN - PIK3CA/PIK3R1	Full-length	1-1068(end)	NP_006209.2	N-terminal DYKDDDDK	128 kDa	Insect (<i>Sf21</i>)
01-438-20N	BTN - PKC η (PRKCH)	Full-length	1-683(end)	NP_006246	N-terminal DYKDDDDK	82 kDa	Insect (<i>Sf21</i>)
01-444-20N	BTN - PKN1	Full-length	1-942(end)	NP_002732.3	N-terminal DYKDDDDK	108 kDa	Insect (<i>Sf21</i>)
05-457-20N	BTN - PLK1	Full-length	1-603(end)	NP_005021.2	N-terminal DYKDDDDK	72 kDa	Insect (<i>Sf21</i>)
08-438-20N	BTN - PYK2(PTK2B)	Full-length	1-967(end)	NP_775267.1	N-terminal DYKDDDDK	115 kDa	Insect (<i>Sf21</i>)
08-459-20N	BTN - RET	Cytoplasmic domain	658-1114(end)	NP_066124.1	N-terminal DYKDDDDK	56 kDa	Insect (<i>Sf21</i>)
01-409-20N	BTN - ROCK1	Catalytic domain	1-477	NP_005397.1	N-terminal DYKDDDDK	59 kDa	Insect (<i>Sf21</i>)
11-405-20N	BTN - SPHK1	Full-length	1-384(end)	NP_001136074.1	N-terminal DYKDDDDK	46 kDa	Insect (<i>Sf21</i>)
08-473-20N	BTN - SRC	Full-length	1-536(end)	NP_005408.1	N-terminal DYKDDDDK	64 kDa	Insect (<i>Sf21</i>)
04-460-20N	BTN - SRPK1	Full-length	1-654(end)	CAI20544.1	N-terminal DYKDDDDK	78 kDa	Insect (<i>Sf21</i>)
11-406-20N	BTN - SRPK2	Full-length	1-618(end)	NP_001191089.1	N-terminal DYKDDDDK	69 kDa	Insect (<i>Sf21</i>)
08-476-20N	BTN - SYK	Full-length	1-635(end)	NP_003168.2	N-terminal DYKDDDDK	76 kDa	Insect (<i>Sf21</i>)
09-419-21C	BTN - TAK1-TAB1(MAP3K7)	Fused protein	1-303	NP_663304.1	C-terminal DYKDDDDK	45 kDa	Insect (<i>Sf21</i>)
09-441-20N	BTN - TGFBR1(TGFBR1)	Catalytic domain	200-503(end)	NP_004603.1	N-terminal DYKDDDDK	38 kDa	Insect (<i>Sf21</i>)
08-485-20N	BTN - TIE2	Cytoplasmic domain	771-1124(end)	NP_000450.1	N-terminal DYKDDDDK	44 kDa	Insect (<i>Sf21</i>)
07-438-20N	BTN - TNIK	Catalytic domain	1-314	NP_055843.1	N-terminal DYKDDDDK	39 kDa	Insect (<i>Sf21</i>)
08-486-20N	BTN - TRKA(NTRK1)	Cytoplasmic domain	436-790(end)	NP_001012331.1	N-terminal DYKDDDDK	44 kDa	Insect (<i>Sf21</i>)
08-487-20N	BTN - TRKB(NTRK2)	Cytoplasmic domain	456-822(end)	NP_001018074.1	N-terminal DYKDDDDK	45 kDa	Insect (<i>Sf21</i>)
08-497-20N	BTN - TRKC(NTRK3)	Catalytic domain	456-825(end)	NP_002521.2	N-terminal DYKDDDDK	46 kDa	Insect (<i>Sf21</i>)
08-447-20N	BTN - TYK2	Catalytic domain	871-1187(end)	NP_003322.3	N-terminal DYKDDDDK	40 kDa	Insect (<i>Sf21</i>)
05-479-20N	BTN - WNK1	Catalytic domain	1-491	NP_061852.1	N-terminal DYKDDDDK	59 kDa	Insect (<i>Sf21</i>)
05-482-20N	BTN - WNK4	Catalytic domain	1-444	NP_115763.2	N-terminal DYKDDDDK	53 kDa	Insect (<i>Sf21</i>)

Updated: 2015/4/14

Note: The contents of this list are subject to change without notice.