



PRODUCTS AND SERVICES

*Drug Discovery Tools
from Carna Biosciences*

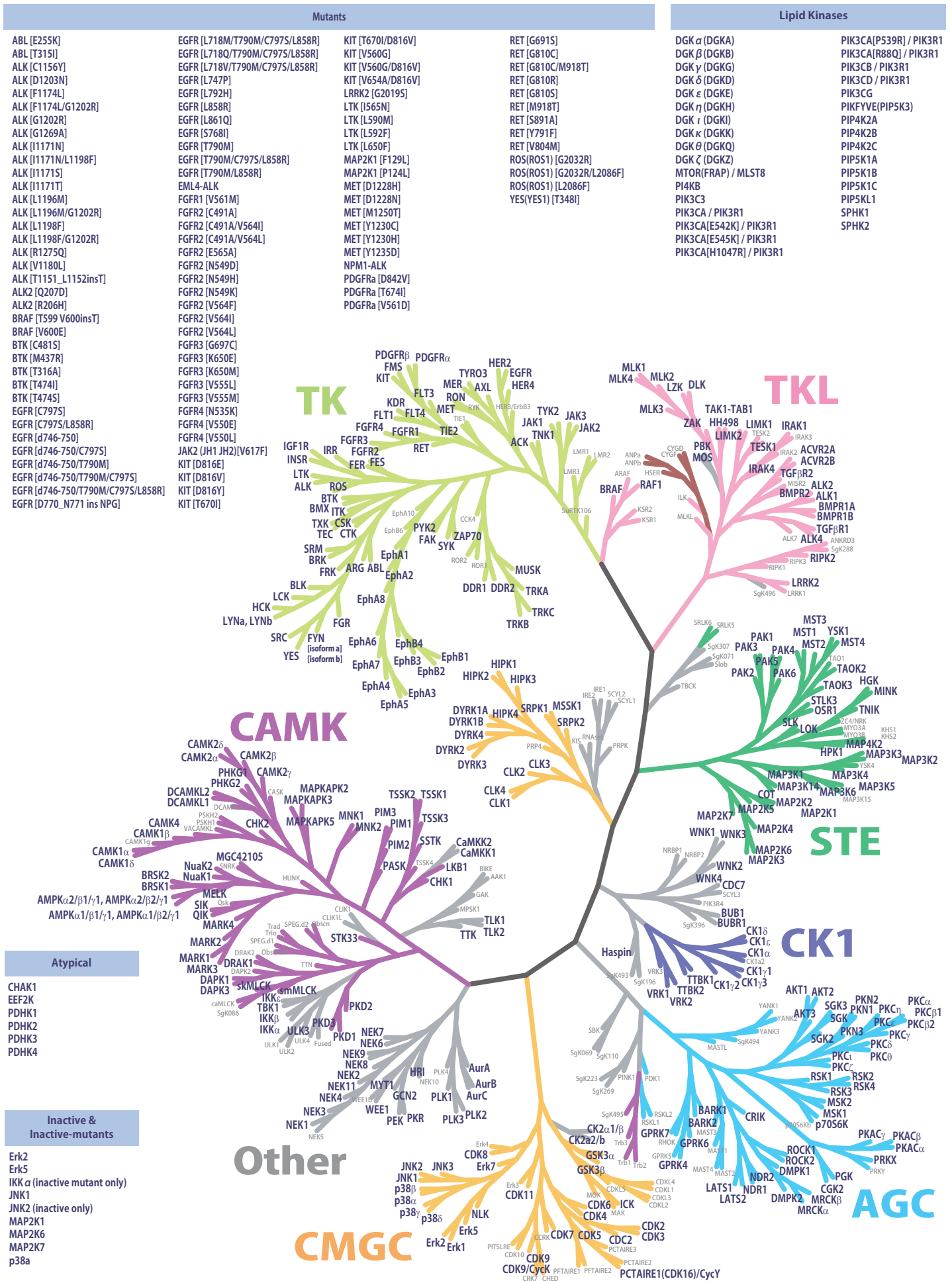
~your kinase company~



Kinases ~Available from Carna Biosciences~	P.1
High Quality Active Kinases	P.2
Biotinylated Kinases	P.3
QuickScout Screening Assist™ Kits	P.4
Biochemical Kinase Screening and Profiling Services	P.5
• Mobility Shift Assay	
• IMAP™ Assay	
• ADP-Glo™ Assay	
• Preincubation Kinase Profiling Service	
• QuickScout™ Pre-selected Panel Profiling Options	
> MAPK Cascade Panel	
> Tyrosine Kinase(TK) Panel	
> Serine/Threonine Kinase (STK) Panel	
> Cell Cycle Panel	
NanoBRET™ TE Intracellular Kinase Cell-Based Assay	P.7
• IC ₅₀ determinations	
• Residence Time Analysis	
• CDK Panel Assay Service	
• Kinome-Wide Profiling Service(K192)	
Detection of Protein-Protein Interactions ~ProbeX™~	P.9
Products & Services by Kinase List	P.10



Kinases ~Available from Carna Biosciences~



High Quality Active Kinases

Carna offers an expanded portfolio of high purity, active human kinases and related products. All of our >450 kinases and products are developed and produced entirely in-house from gene cloning, expression and purification, and undergo rigorous quality control. All products are available in various sizes starting from 5µg up to bulk production, are delivered to you with a lot specific data sheet.

*Target lists: Please refer to the attachment.

Industry Leading Production and Quality Control Methods

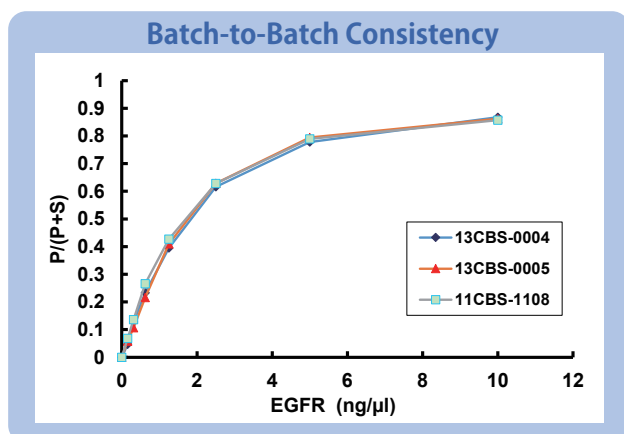
- Proprietary Production Methods ensure batch-to-batch consistency (Fig.1)
- Relevant kinase specific substrate(s) measure activity of each batch
- Target kinase activity maximized to minimize literature documented host cell activity
- Kinase constructs are carefully selected from published literature
- DNA sequence confirmed prior to protein expression
- Amino acid sequence confirmed by Peptide Mass Fingerprinting(PMF)

Unparalleled Kinase Activity

Production process yields highly active kinases.

Some kinases are activated further by :

- ◆ Expression with upstream kinase(s)
- ◆ Tag removal
- ◆ ATP treatment



[Fig.1]

Lipid Kinases and Related Products

Carna's lipid kinase panel includes numerous members of the PIK family. The enzymes enable complete and easy investigation of lipid kinase drug targets.

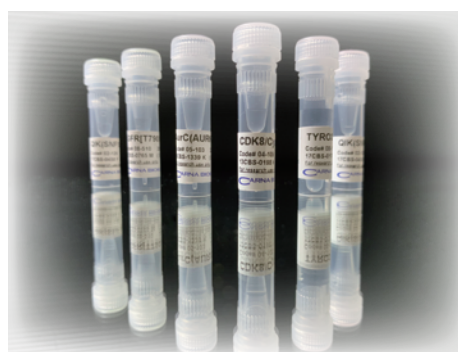
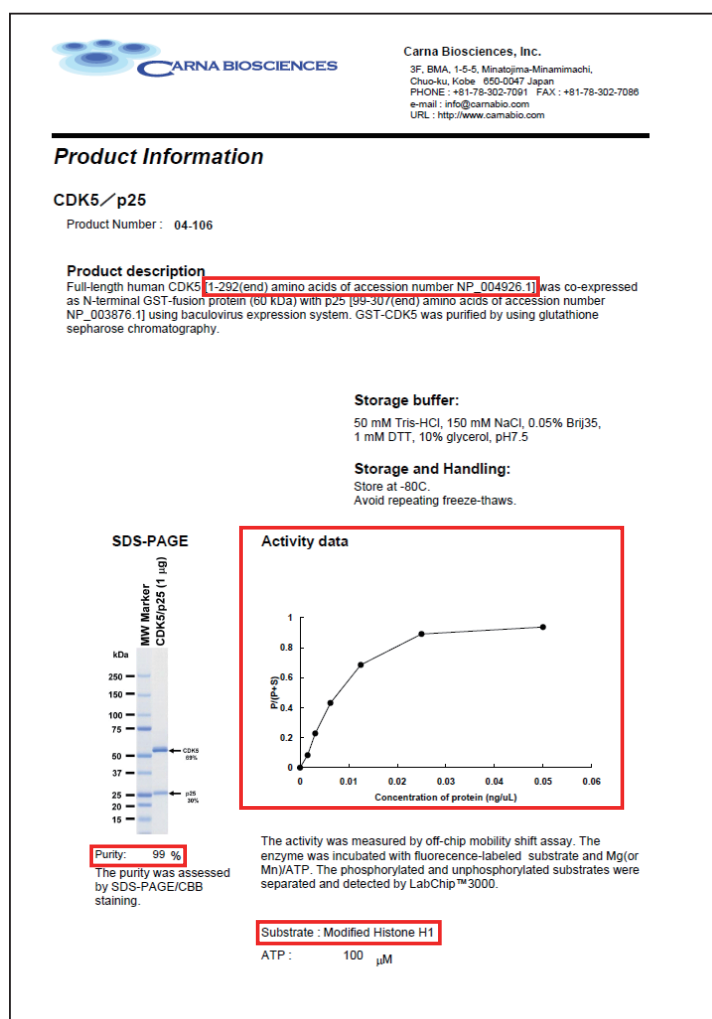
Protein Substrates

Our protein substrates are available for use in phosphorylation activity assays to assess kinase activity.

Custom protein production

We offer custom production services using our extensive expertise, for any kinase of interest.

For more information, please contact us at info@carnabio.com



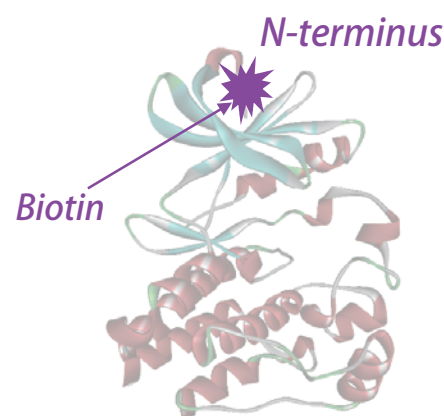
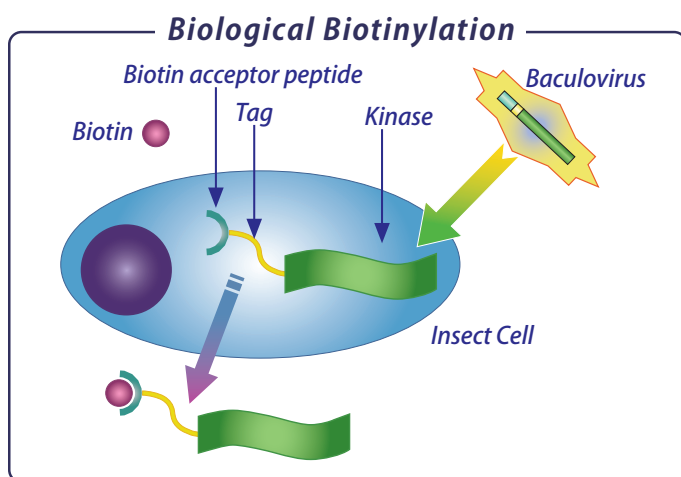
Biotinylated Kinases

Biotinylated Kinases are best suited for the study of compound binding affinity and other kinase-molecule interactions using devices measuring Surface Plasmon Resonance (SPR), BioLayer Interferometry (BLI) and other similar biomolecular interactions. They can also be utilized in homogenous proximity-based binding assays such as TR-FRET, AlphaScreen™ and HTRF® to interrogate inhibitor binding affinity, determine on-off rates, and measure binding kinetics. The immobilization of target proteins onto sensor surfaces without impairing their structure and activity can be challenging in small molecule drug discovery. Carna's in-house, single-site specifically biotinylated kinases are easily immobilized, leading to rapid acquisition of accurate and real-time data for evaluation of your drug candidates!

*Target lists: Please refer to the attachment.

Advantages of Carna's Biological Biotinylation Process

- Kinases are labeled with a single biotin at the N-terminus
- Easy-to-use ; no additional labeling required
- Native, catalytically active kinase domains are preserved
- High quality human kinases produced via Baculovirus expression system
- Stable activity determined post expression
- Select kinases available pre-activated (via ATP treatment) and non-activated (without ATP treatment)



SPR data samples* using Carna's biotinylated kinases are shown on our website

(*Measured using Biacore T200 in collaboration with Oncolines B.V)

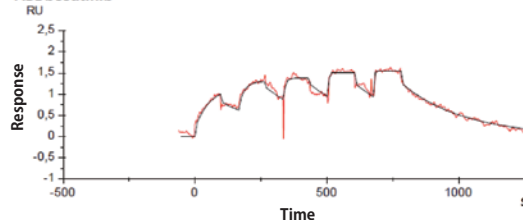
Biotinylated Kinases				
Catalog No.	Product Name	Product Size		
08-401-20N	BTN-ABL(ABL1)	10ug	100ug	Bulk
01-401-20N	BTN-AKT1	10ug	100ug	Bulk
01-402-20N	BTN-AKT2	10ug	100ug	Bulk
08-405-20N	BTN-ALK	10ug	100ug	Bulk
08-429-20N	BTN-ALK[L1196M]	10ug	100ug	Bulk
05-401-20N	BTN-AurA (AURKA)	10ug	100ug	Bulk
05-402-21N	BTN-AurB (AURKB)/INCENP*	10ug	100ug	Bulk
08-407-20N	BTN-AXL	10ug	100ug	Bulk
08-479-20N	BTN-BMX	10ug	100ug	Bulk
09-422-20N	BTN-BRAF	10ug	100ug	Bulk
08-480-20N	BTN-BTK *	10ug	100ug	Bulk
08-417-20N	BTN-BTK[C481S] *	10ug	100ug	Bulk
08-417-23N	BTN-BTK[C481S][non-activated]	10ug	100ug	Bulk
08-480-23N	BTN-BTK[non-activated]	10ug	100ug	TBA
08-418-20N	BTN-BTK[T316A] *	10ug	100ug	Bulk
08-418-23N	BTN-BTK[T316A][non-activated]	10ug	100ug	Bulk
08-419-20N	BTN-BTK[T474I] *	10ug	100ug	Bulk
08-419-23N	BTN-BTK[T474I][non-activated]	10ug	100ug	Bulk
08-420-20N	BTN-BTK[T474S] *	10ug	100ug	Bulk

Biacore Sensorgram
(Reference Data)

Click!

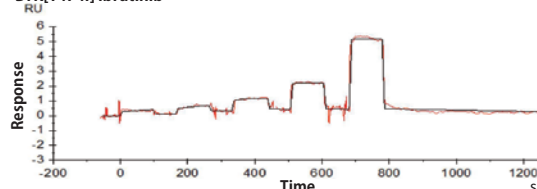
Binding kinetics of bosutinib to ABL

ABL bosutinib



Binding kinetics of ibrutinib to BTK[T474I]

BTK[T474I] Ibrutinib



QuickScout Screening Assist™ Kits

QuickScout Screening Assist™ Kits are designed to accelerate your in house compound screening, particularly secondary and counter- screening applications. Our Kits provide essential reagents and detailed assay protocols in one package, and are available for more than 300 human kinase targets. After your initial kit purchase, components contained in the kit can be purchased separately and in bulk.

*Target lists: Please refer to the attachment.

Advantages of Carna's Assay Kits

- Prepared utilizing the extensive expertise of our kinase profiling team
- Ready-To-Run products & protocols save time and money
- Scalable for HTS applications

Designed for primary, in-house screening procedures applicable to Lead Generation through Lead Optimization!



Assay Platform	Minimum Kit Size	Kit Components
Mobility Shift Assay QSS Assist™ MSA This MSA kit works best using LabChip® technology from PerkinElmer, Inc.	400dp Equivalent to 1 x 384-well plate	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture (ATP, Cation included) ● Assay Buffer ● Termination Buffer ● Assay Protocol (Separation conditions included)
FP(IMPACT™) QSS Assist™ FP	400dp Equivalent to 1 x 384-well plate	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture (ATP, Cation included) ● Assay Buffer ● Assay Protocol
TR-FRET QSS Assist™ TR-FRET	400dp Equivalent to 1 x 384-well plate	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture (ATP, Cation included) ● Assay Buffer ● Assay Protocol
ELISA QSS Assist™ ELISA	100dp Equivalent to 1 x 96-well plate	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture (ATP, Cation included) ● Assay Buffer ● Antibody for ELISA (except for TTK & WEE1) ● Assay Protocol
ADP-Glo™ QSS Assist™ ADP-Glo™	400dp Equivalent to 1 x 384-well plate	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Solution ● Kinase dilution Buffer (Cation included) ● Assay Buffer ● MgCl2 solution for detection reagent ● Assay Protocol

Sample protocols for all kit platforms are available online.
 Each kinase kit is made-to-order, with turnaround time of 2-3 weeks.

Biochemical Kinase Screening and Profiling Services

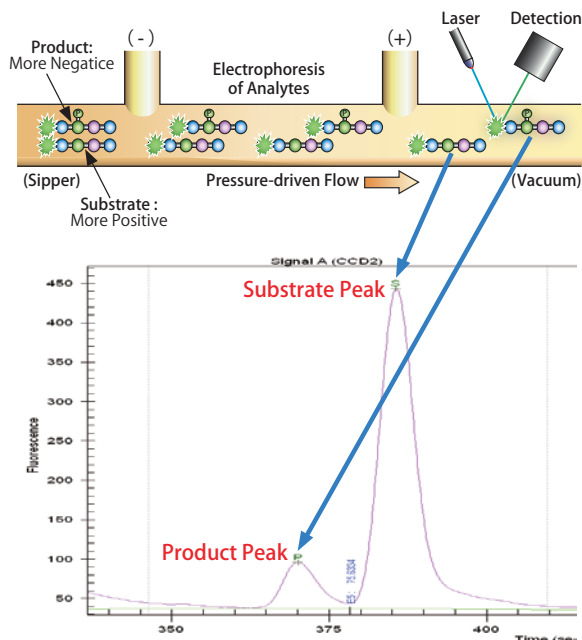
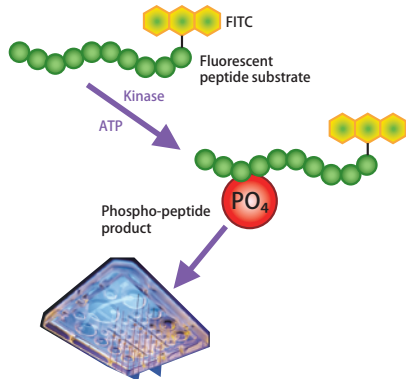
Carna offers >320 biochemical screening and profiling assays for assessing potency and selectivity of your compounds. Our three assay platforms allow us to interrogate a wide range of kinase targets using stringent SOPs. We offer testing at higher ATP concentration (1mM) to provide insight into compound inhibition and potency under more physiologically relevant conditions, in addition to our ~ATP km assay services.

*Target lists: Please refer to the attachment.

Three Assay Platforms

1. Mobility Shift Assay

Direct monitoring of phosphorylation by measuring non-phosphorylated and phosphorylated substrate.



2. IMATM (Immobilized Metal Ion-Affinity Partitioning) Assay

3. ADP-GloTM Assay

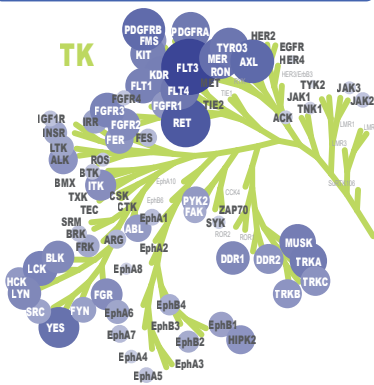
~ATP Km / 1mM ATP Assays

ATP concentration approximating Km are our routine kinase assays

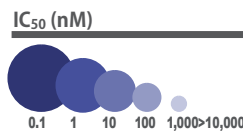
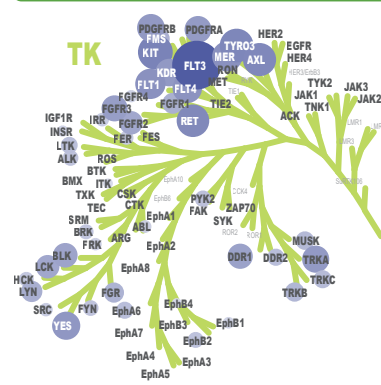
Many assays are also available at 1mM ATP

Assays approaching physiological ATP levels, provide insight to in vivo pharmacology.

Sunitinib vs. TKs at [ATP]=Km bin.

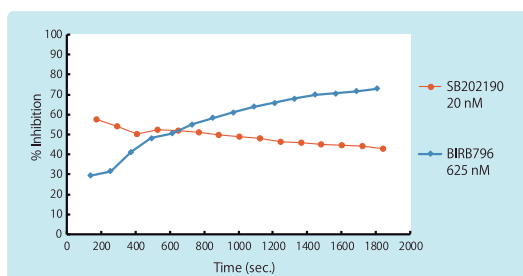


Sunitinib vs. TKs at [ATP]=1mM



Preincubation Kinase Profiling Service

Carna's preincubation service can be utilized to study slow binding compounds. This service incorporates a thirty (30) minute room temperature preincubation of target kinase with your test compound(s) prior to measuring activity in our standard Mobility Shift Assay.



Time course of p38α inhibitors (ATP=1mM)

Structurally unrelated SB202190 and BIRB796 are potent inhibitors of p38α. BIRB796 interacts with p38α in a manner different from SB202190, and its binding induces a slow conformational change that locks the protein into an inactive conformation. The potency of BIRB796 increases with prolonged incubation, which is easily detected utilizing the pre-incubation service.

Biochemical Kinase Screening and Profiling Services

QuickScout™ Pre-selected Panel Profiling Options

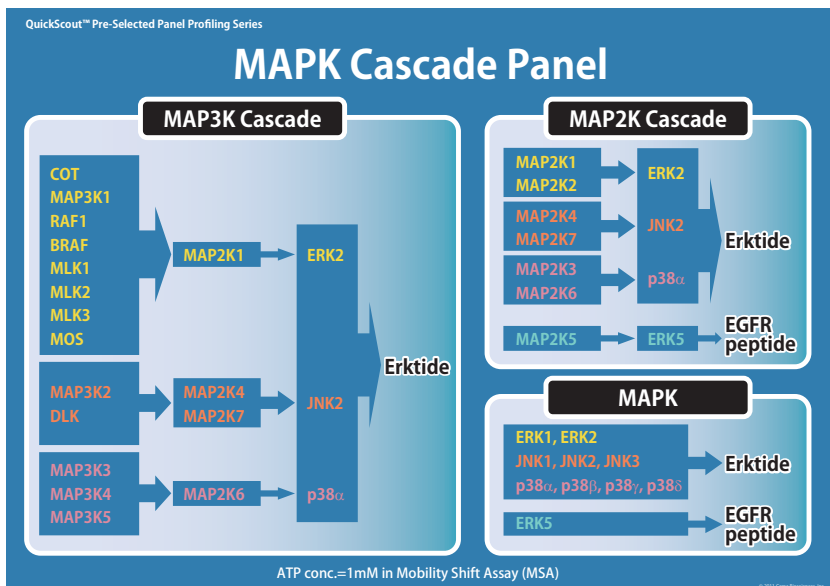
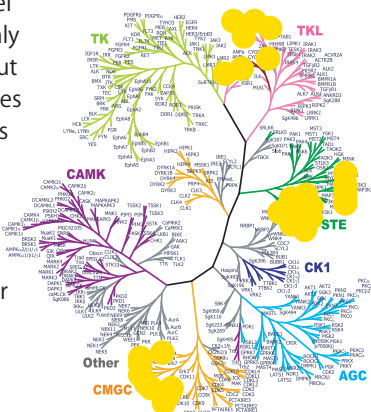
Our selected panel series are well-suited for an initial profiling of your compounds.

*Target lists: Please refer to the attachment.

MAPK Cascade Panel ver. 2.0

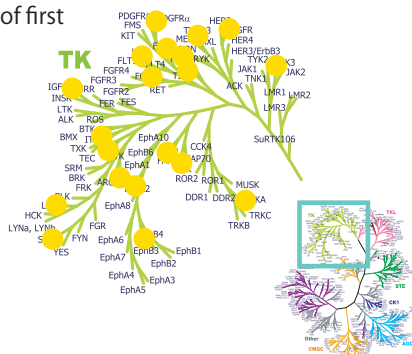
The MAPK cascade plays an important role in the intracellular signal transduction of eukaryotic cells.

Our MAPK panel includes not only MAP kinases, but upstream kinases such as MAPKKs and MAPKKKs. This panel is useful for analyzing the function of your compounds in the cascade.



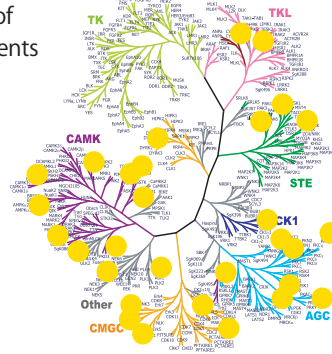
Tyrosine Kinase (TK) Panel ver. 2.0

QuickScout™ TK Panel consists of 20 Pre-Selected receptor and non-receptor Tyrosine Kinases, and helps you to rapidly screen your compounds against druggable and clinically relevant kinases. Identification and optimization of small molecule inhibitors against many of the targets in our 20 TK Panel have led to the development of first and second generation therapeutics for the treatment of diseases such as leukemia and cancers of the lung, breast, and kidney.



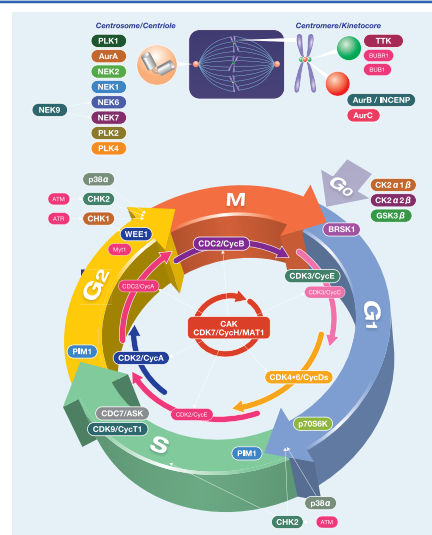
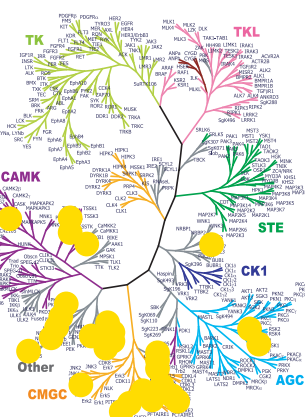
Serine/Threonine Kinase (STK) Panel ver. 3.0

QuickScout™ STK Panel consists of 30 Pre-Selected Serine / Threonine kinases that are key members of the AGC, CAMK, CMGC, STE, TKL, and Other Group of STK kinases. This Panel allows you to screen your lead compounds using the industry's most diverse Kinome Sampler and helps to discover and characterize the selectivity of compounds as potential treatments for cancer, inflammatory, metabolic and/or neurological diseases.



Cell Cycle Panel ver. 2.0

QuickScout™ Cell Cycle Panel is comprised of relevant kinases for cell-cycle regulation and is well-suited for determining whether your compound acts on cell division. This panel mainly includes kinases that are directly involved in the cell-cycle where their inhibition may interfere with cell proliferation.



NanoBRET™ TE Intracellular Kinase Cell-Based Assay

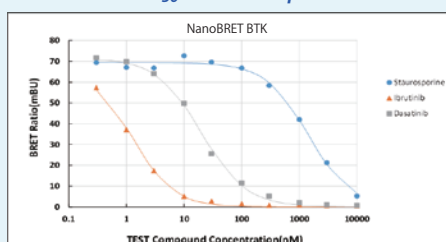
Quantifying kinase inhibitor occupancy, selectivity, and affinity within the cellular environment is crucial to more accurately predict engagement potencies against human kinases. In addition to equilibrium evaluation, kinetic parameters such as Residence Time should be determined for better compound optimization. Quantitative and wide-spectrum kinase profiling services using the NanoBRET™ Target Engagement Intracellular Kinase Assay System (Promega) enable you to assess your compound's engagement for a selected intracellular target under physiological conditions, including compound Residence Time at the target, while keeping the cells intact. Simply submit your compound(s) of interest, and Carina will rapidly deliver cellular IC₅₀ values and Residence Time!

*Target lists: Please refer to the attachment. For targets not listed, please inquire.

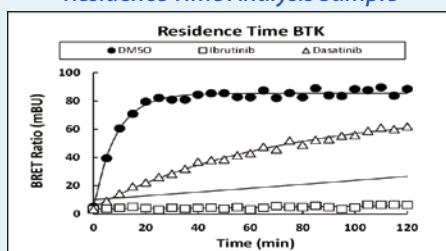


● Investigate Inhibitor Binding and Residence Time in Intact Cells Expressing Full-length Kinases

IC₅₀ Data Sample



Residence Time Analysis Sample



IC₅₀ Determinations & Residence Time Analysis

IC₅₀ determinations

(7 serial half-log dilutions, 8 conc. points)

Turnaround:
2 weeks upon receipt of compounds

Residence Time Analysis*

(performed in duplicate based on your IC₅₀ evaluation result)

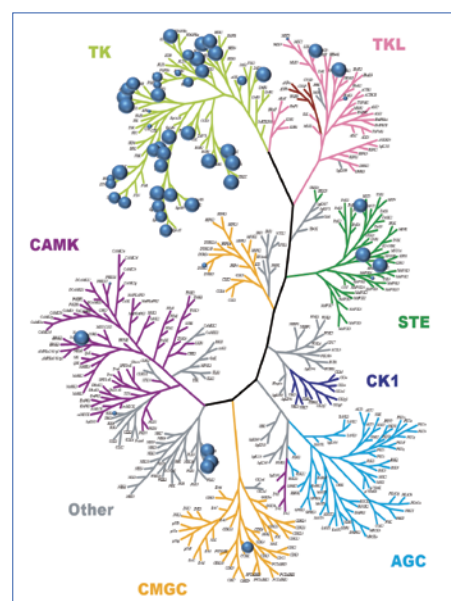
Turnaround: 3 weeks after dose setting
*Preliminary IC₅₀ determinations required for study dose setting.

Panel Services

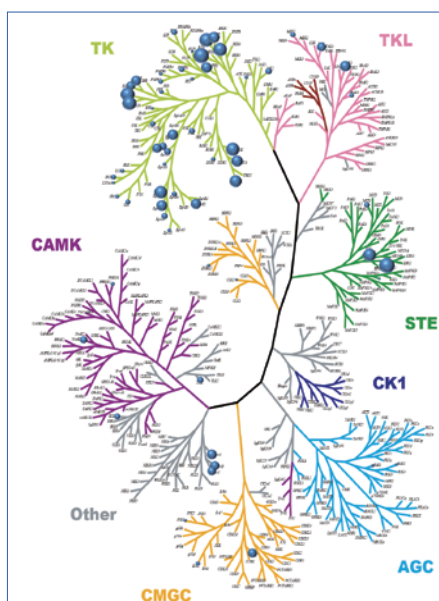
- ◆ CDK Panel Assay Service
- ◆ Kinome-Wide Profiling Service (192 Kinase Panel)

● Compare and Assess Selectivity and Potency at Cellular ATP Concentration

[Target occupancy using 1 μ M Crizotinib]



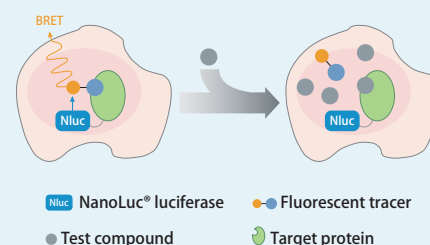
Carina Kinase Profiling service
ATP Km, Mobility Shift Assay



NanoBRET™ (HEK 293 cells)
Cell Chem Biol. 2018 Feb 15;25(2):206-214.e11

NanoBRET™ System

A cell-permeable fluorescent NanoBRET™ tracer, a BRET acceptor, is added to HEK293 cells expressing a full length kinase/Nano-Luc® fusion protein. Engagement of the tracer to the target protein generates a BRET signal.



Binding of the test compound to the target protein results in a loss of NanoBRET™ signal between the target protein and the tracer inside intact cells.

NanoBRET™ TE Intracellular Kinase Cell-Based Assay

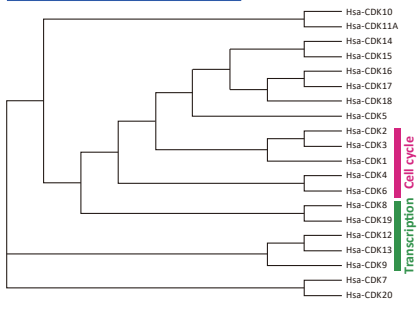
Two Panel Assay Services using NanoBRET™ Target Engagement (TE) Intracellular Kinase Assay technology are available to accelerate your investigations.

*Target lists: Please refer to the attachment. For targets not listed, please inquire.

● CDK Panel Assay Service

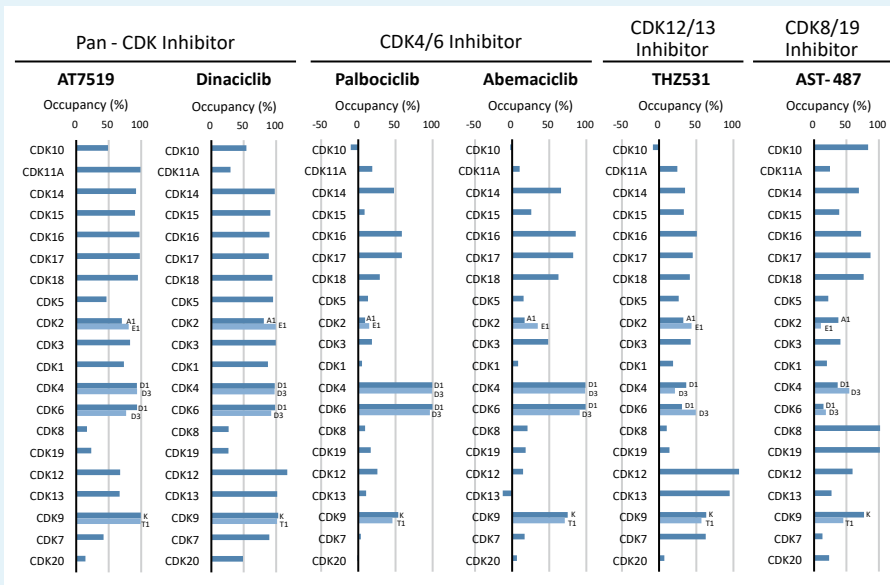
Carna's NanoBRET™ TE Intracellular CDK Panel Assay consists of 24 CDK targets. It is an ideal tool to examine the specificity of your CDK inhibitor across an extensive set of CDK family kinases: in the presence of the cyclin subunit; in a cellular environment; and in a unified assay format. The compound's engagement with each CDK is quantitatively measured inside living HEK293 cells, which provides a snapshot of your compound's selectivity across all the CDK targets in the panel.

CDK Phylogenic Tree



(Right)

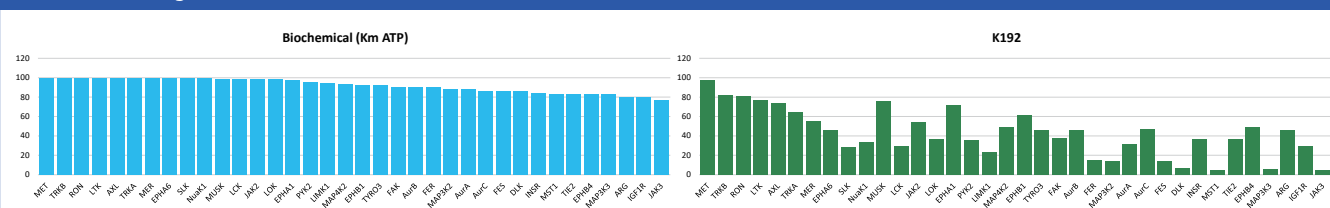
NanoBRET™ TE Intracellular CDK Panel Assay reveals each selectivity at 10 μ M conc. of pan-CDK inhibitors (AT7519, dinaciclib), CDK4/6 inhibitors (palbociclib, abemaciclib), CDK12/13 inhibitor (THZ531), and CDK8/19 inhibitor (AST-487).



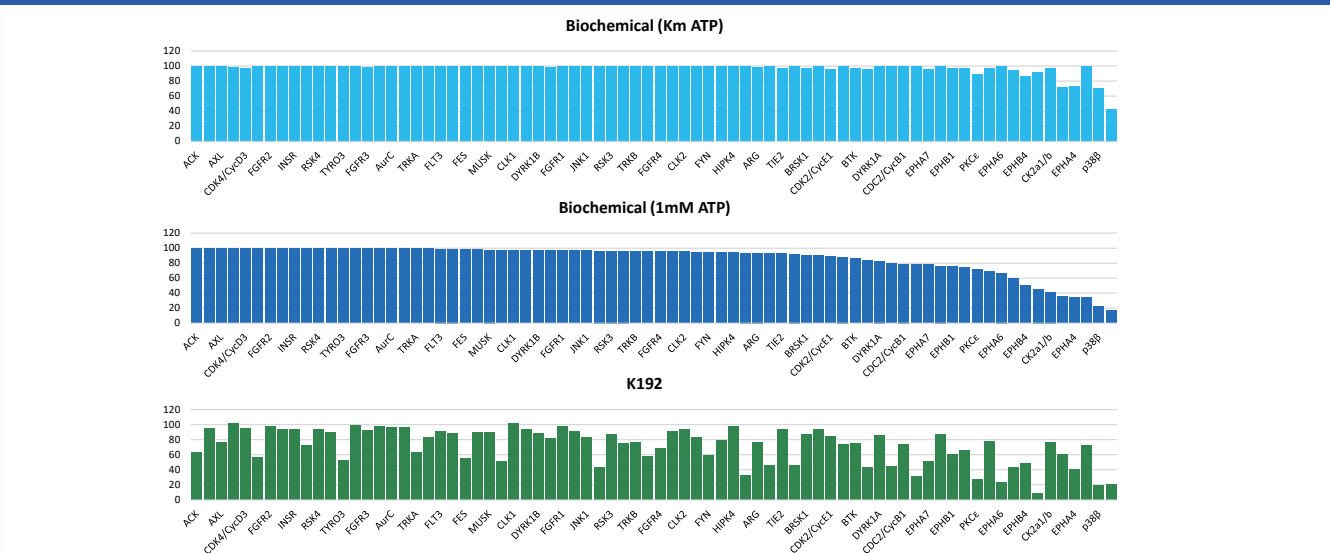
● Kinome-Wide Profiling Service (K192)

NanoBRET™ TE Intracellular Kinase Assay System (Promega) is used in this broad panel of 192 kinase targets (see list), each expressed intracellularly by transient transfection. Compound evaluation against all targets is performed simultaneously at 1 compound concentration in duplicate, under the same assay conditions. Follow up IC₅₀ determinations also available.

Comparison data using 1 μ M Crizotinib in a biochemical assay (Mobility Shift Assay) and NanoBRET™ K192 cell assay panel (36 selected targets shown)



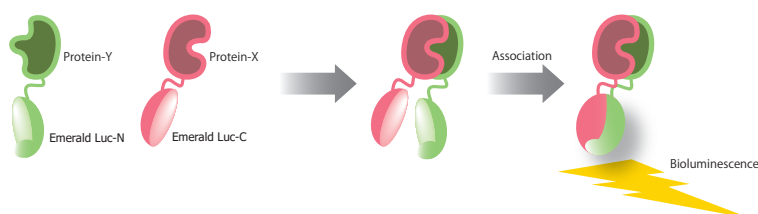
Comparison data using 0.3 μ M CC1 inhibitor in a biochemical assay (Mobility Shift Assay) and NanoBRET™ K192 cell assay panel



Detection of Protein-Protein Interactions ~ProbeX™~

Carna's split luciferase complementation assay, utilizing a unique luciferase derived from *Pyrearinus termitilluminans* (Emerald Luciferase, E-Luc), is a valuable tool for your study of Protein-Protein Interactions (PPIs). Detection of various types of PPIs, including GPCRs, is performed with ease and high-sensitivity. In addition to off-the-shelf cell lines, we develop stable transfected cell lines suitable for detecting specific PPIs of interest.

● Split Luciferase Technology



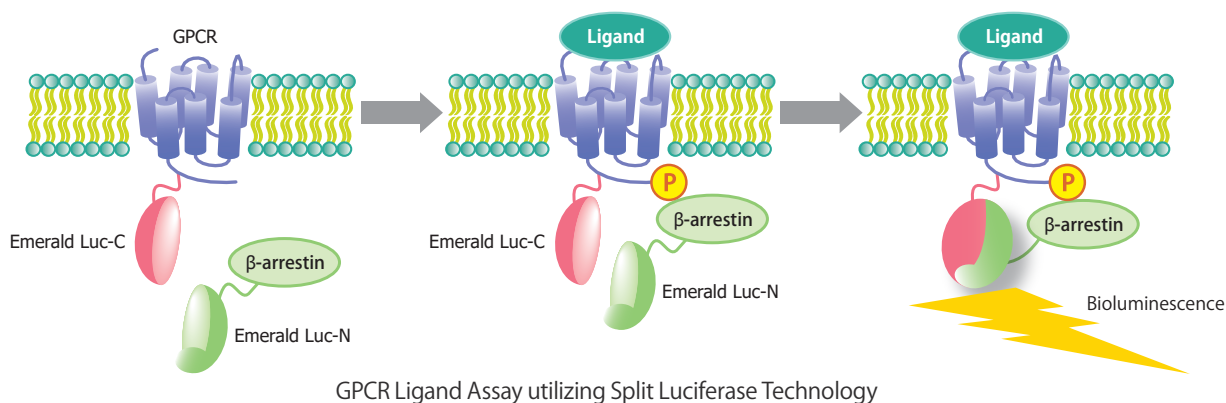
E-Luc is known to emit a brighter and more stable signal than conventional firefly luciferases. The N- and C-terminal domains of luciferase can be separated into two fragments, which can then re-associate in cells. When the two fragments of the reporter proteins are brought within proximity, they spontaneously refold and generate a detectable signal (patent filed).

● Application for GPCR

The N-terminal and C-terminal fragments of emerald split luciferase are fused to β -arrestin and GPCR, respectively. Binding of a ligand to the GPCR triggers phosphorylation of the GPCR, thereby inducing its interaction with β -arrestin. This interaction brings the N-terminal luciferase fragment into proximity with the C-terminal fragment, and bioluminescence activity is recovered.

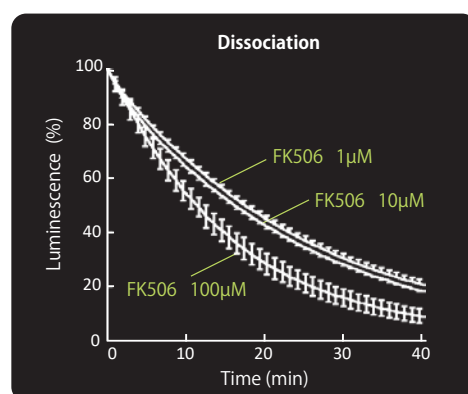
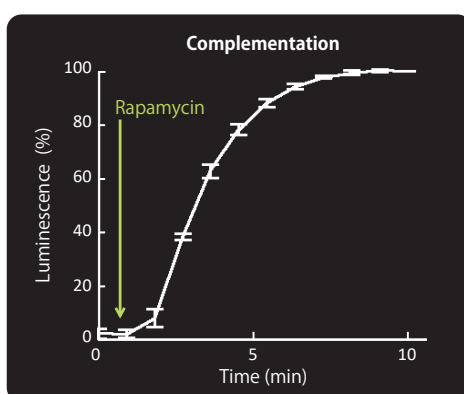
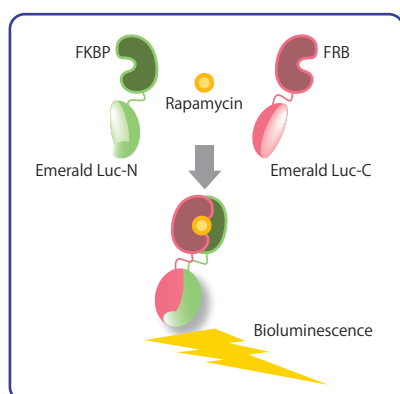
Please view the list* showing our validated stable transfectants. We also develop custom cell lines suitable for your needs on a fee for service basis.

*Target lists: Please refer to the attachment.



● Example Application - FKBP-FRB Interaction

The immunosuppressant macrolide, rapamycin, mediates the interaction of two proteins: FKBP and FRB (left). Addition of rapamycin facilitates E-Luc complementation and a concomitant rapid increase in bioluminescence (middle). A competitive inhibitor of FK506 decoupled the rapamycin-induced signal (right), indicating the E-Luc complementation is reversible. This is in contrast to the split GFP system which is irreversible and not applicable to detect dissociation of two proteins.



		Products						Services								
								Cell-Free Assays				Cell-Based Assays				
<div>Products & Services</div>	Kinase Proteins	Biotiny-lated Kinases	Kinase Protein Assay Kit					Mobility Shift Assay/IMAP™				ADP-Glo™ (ATP=Km)	NanoBRET™ TE Intracellular Kinase Cell-Based Assay Services			
								MSA/IMAP™		ATP conc.						
			Kinases name			MSA	FP (IMAP™)	ELISA	TR-FRET	ADP-Glo™	■: IMAP™ (FP)		Preincubation Study	Km	1mM	
AAK1														○		○
ABL(ABL1)	○	○	○						○	○	○	○		○		
ABL(ABL1) [E255K]	○		○						○	○	○	○		○		
ABL(ABL1) [F317I]														○		
ABL(ABL1) [F317L]														○		
ABL(ABL1) [H396P]														○		
ABL(ABL1) [M351T]														○		
ABL(ABL1) [Q252H]														○		
ABL(ABL1) [T315I]	○	○	○						○	○	○	○		○		
ABL(ABL1) [Y253F]														○		
ACK(TNK2)	○		○						○		○	○		○		○
ACTR2B(ACVR2B)	○	○				○										
ACVR1C(ALK7)	○															
ACVR2A	○	○				○										
ACVRL1(ALK1)	○					○								○		
ADK														○		
AKT1	○	○	○	○					○	○	○	○		○		
AKT1 [E17K]														○		
AKT2	○	○	○	○					○	○	○			○		○
AKT2 [E17K]														○		
AKT3	○		○	○					○	○	○					
AKT3 [E17K]														○		
AKT3 [G171R]														○		
ALK	○	○	○			○			○	○	○	○		○		
ALK [C1156Y]	○	○	○						○	○	○	○				
ALK [D1203N]	○															
ALK [F1174L]	○	○	○						○	○	○	○				
ALK [F1174L/G1202R]	○	○														
ALK [G1202R]	○	○	○						○	○	○	○				
ALK [G1269A]	○	○	○						○	○	○	○				
ALK [I1171N]	○	○														
ALK [I1171N/L1198F]	○	○														
ALK [I1171S]	○	○														
ALK [I1171T]	○	○														
ALK [L1196M]	○	○	○						○	○	○	○				
ALK [L1196M/G1202R]	○	○														
ALK [L1198F]	○	○														
ALK [L1198F/G1202R]	○															
ALK [R1275Q]	○	○	○						○	○	○	○				
ALK [T1151_L1152insT]	○		○						○		○	○				
ALK [V1180L]	○	○														
ALK2(ACVR1)	○					○								○		
ALK2(ACVR1) [G328V]																

Products & Services by Kinase

<div>Products & Services</div> <div>Kinases name</div>	Products							Services						
	Kinase Proteins	Biotiny-lated Kinases	Kinase Protein Assay Kit					Cell-Free Assays				Cell-Based Assays		
								Mobility Shift Assay/IMAP™		ADP-Glo™ (ATP=Km)	NanoBRET™ TE Intracellular Kinase Cell-Based Assay Services	Assay Service	CDK Panel	K192 Panel
								MSA/IMAP™	ATP conc.					
			MSA	FP (IMAP™)	ELISA	TR-FRET	ADP-Glo™	■: IMAP™ (FP)	Preincubation Study	Km	1mM			
BTK [C481S]	○	○	○					○		○	○	○		
BTK [C481S] [non-activated]		○												
BTK [E41K]												○		
BTK [M437R]	○	○												
BTK [M437R] [non-activated]		○												
BTK [P190K]												○		
BTK [T316A]		○												
BTK [T316A] [non-activated]		○												
BTK [T474I]		○										○		
BTK [T474I] [non-activated]		○												
BTK [T474S]		○												
BTK [T474S] [non-activated]		○												
BUB1/BUB3	○	○	○					○	○	○				
BUBR1(BUB1B)	○													
CaMK1α(CAMK1)	○		○					○		○	○	○		○
CaMK1β(PNCK)	○											○		
CaMK1γ(CAMK1G)												○		
CaMK1δ(CAMK1D)	○		○					○		○		○		
CaMK2α(CAMK2A)	○		○	○				○	○	○	○	○		○
CaMK2β(CAMK2B)	○	○	○					○	○	○	○			
CaMK2γ(CAMK2G)	○		○					○	○	○	○	○		
CaMK2δ(CAMK2D)	○	○	○					○	○	○	○	○		○
CaMK4	○		○	○				○	○	○	○			
CAMKK1	○													
CAMKK1(124-411)												○		
CAMKK2	○													
CAMKK2(165-445)												○		
CDC7/ASK	○		○					○		○	○			
CDK1(CDC2)/CycA2	○	○										○		
CDK1(CDC2)/CycB1	○	○	○	○				○	○	○	○	○	○	○
CDK1(CDC2)/CycE1												○		
CDK2/CycA1												○	○	
CDK2/CycA2	○	○	○	○				○	○	○	○	○		
CDK2/CycE1	○	○	○					○	○	○	○	○	○	○
CDK3/CycE1	○	○	○	○				○	○	○	○	○	○	○
CDK4/CycD1												○	○	
CDK4/CycD3	○	○	○	○				○		○	○	○	○	○
CDK5												○		
CDK5/p25	○	○	○	○				○	○	○	○			
CDK5/CDK5R1												○	○	○
CDK5/CDK5R2												○		
CDK6/CycD1	○											○	○	○
CDK6/CycD3	○		○					○		○	○	○	○	
CDK7												○		○
CDK7/CycH												○	○	
CDK7/CycH/MAT1	○	○	○					○		○	○			
CDK8/CycC	○	○			○							○	○	
CDK9/CycK	○	○										○	○	○
CDK9/CycT1	○	○	○					○		○	○	○	○	
CDK9/CycT2	○	○												
CDK10/Cycl2												○	○	○
CDK11A/CycK												○		
CDK11A/Cycl2												○	○	
CDK12(CRKR5)/CycK												○	○	
CDK12(CRKR5)(720-1490aa)/CycK		○												
CDK13(CHED)/CycK	○	○										○	○	
CDK13(CHED)(694-1512aa)/CycK		○												
CDK14/CycY												○	○	○
CDK15/CycY												○	○	○
CDK17/CycY												○	○	○
CDK18/CycY												○	○	○
CDK19(CDC2L6)/CycC	○	○			○							○	○	
CDK20/CycH												○	○	○
CDKL1												○		○
CDKL2												○		○
CDKL3												○		○
CDKL5												○		○
CGK2(PRKG2)	○		○	○				○	○	○		○		
CHAK1(TRPM7)	○													
CHK1(CHEK1)	○	○	○	○				○	○	○	○	○		
CHK2(CHEK2)	○		○	○				○	○	○	○	○		○
CK1α(CSNK1A1)	○		○					○		○	○			
CK1α1L(CSNK1A1L)												○		○
CK1γ1(CSNK1G1)	○		○					○	○	○				
CK1γ2(CSNK1G2)	○		○					○	○	○		○		○
CK1γ3(CSNK1G3)	○		○					○	○	○				
CK1δ(CSNK1D)	○		○	○				○	○	○	○	○		○

Products & Services by Kinase

	Products							Services							
<div>Products & Services</div>	Kinase Proteins	Biotiny-lated Kinases	Kinase Protein Assay Kit					Cell-Free Assays				Cell-Based Assays			
								Mobility Shift Assay/ IMAP™			ADP-Glo™ (ATP=Km)	NanoBRET™ TE Intracellular Kinase Cell-Based Assay Services			
			MSA/IMAP™	ATP conc.											
Kinases name			MSA	FP (IMAP™)	ELISA	TR-FRET	ADP-Glo™	■: IMAP™ (FP)	Preincubation Study	Km	1mM		Assay Service	CDK Panel	K192 Panel
CK1ε(CSNK1E)	○		○					○		○	○		○		
CK2α1(CSNK2A1)													○		○
CK2α1/β(CSNK2A1/B)	○		○					○	○	○	○				
CK2α2(CSNK2A2)													○		○
CK2α2/β(CSNK2A2/B)	○		○					○	○	○					
CLK1	○		○	○				○	○	○	○		○		○
CLK2	○		○					○	○	○	○		○		○
CLK3	○		○					○	○	○					
CLK4	○												○		○
Cofilin2	○														
COQ8B													○		
COT(MAP3K8)	○				○			○			○				
CRIK(CIT)	○		○	○				○		○					
CSK	○		○			○		○		○	○		○		
DAPK1	○		○	○				○		○	○				
DAPK2													○		○
DAPK3	○	○													
DCAMKL1(DCLK1)	○												○		
DCAMKL2	○		○					○		○					
DCLK3													○		○
DDR1	○	○	○					○		○	○		○		
DDR1(SRC treated)	○	○													
DDR2	○	○	○					○		○	○		○		
DDR2 [N456S]													○		
DGKa(DGKA)	○	○					○					○			
DGKa(DGKA)(196-735aa)		○													
DGKβ(DGKB)	○	○					○					○			
DGKβ(DGKB)(235-803aa)		○													
mouse DGKβ(DGKB)		○													
DGKy(DGKG)	○	○					○					○			
DGKy(DGKG)(261-791aa)		○													
mouse DGKy(DGKG)		○													
DGKδ(DGKD)	○	○					○					○			
DGKe(DGKE)	○	○					○					○			
DGKη(DGKH)	○	○					○					○			
DGKi(DGKI)	○	○					○					○			
DGKκ(DGKK)	○	○										○			
DGKθ(DGKQ)	○	○					○					○			
DGKζ(DGKZ)	○	○					○					○			
DLK(MAP3K12)	○	○			○			○			○		○		○
DMPK1(DMPK)	○														
DMPK2(CDC42BPG)	○														
DRAK1(STK17A)	○														
DRAK2(STK17B)													○		○
DYRK1A	○		○					○	○	○	○		○		○
DYRK1B	○		○					○	○	○	○		○		○
DYRK2	○		○					○	○	○	○		○		
DYRK3	○		○					○	○	○	○				
DYRK4	○		○												
EEF2K	○		○					○		○					
EGFR(ERBB1)	○	○	○					○	○	○	○		○		
EGFR [C797S]	○	○													
EGFR [C797S/L858R]	○	○						○		○	○				
EGFR [d746-750]	○	○	○					○	○	○	○		○		
EGFR [d746-750/C797S]	○	○						○	○	○	○				
EGFR [d746-750/T790M]	○		○					○		○	○				
EGFR [d746-750/T790M/C797S]	○							○			○				
EGFR [d746-750/T790M/C797S/L858R]	○	○													
EGFR [D770_N771insNPG]	○		○					○		○	○				
EGFR [D770_N771insNPG/T790M]	○														
EGFR [L747P]	○														
EGFR [L792H]	○														
EGFR [L858R]	○	○	○					○	○	○	○		○		
EGFR [L861Q]	○		○					○	○	○	○				
EGFR [S768I]	○														
EGFR [T790M]	○	○	○					○	○	○	○				
EGFR [T790M/L858R]	○	○	○					○	○	○	○		○		
EGFR [T790M/C797S/L858R]	○							○		○	○				
EIF2S1	○														
EIF4EBP1	○														
EML4-ALK	○		○					○		○	○				
EPHA1	○		○					○	○	○	○		○		○
EPHA2	○	○	○			○		○	○	○	○		○		
EPHA2 [non-activated]		○													
EPHA3	○		○					○	○	○	○		○		
EPHA4	○		○					○	○	○	○		○		○
EPHA5	○		○					○	○	○	○		○		

Products & Services by Kinase

Products & Services	Products							Services						
	Kinase Proteins	Biotinylated Kinases	Kinase Protein Assay Kit					Cell-Free Assays				Cell-Based Assays		
								Mobility Shift Assay/IMAP™		ADP-Glo™ (ATP=Km)		NanoBRET™ TE Intracellular Kinase Cell-Based Assay Services		
								MSA/IMAP™	ATP conc.					
			MSA	FP (IMAP™)	ELISA	TR-FRET	ADP-Glo™	■: IMAP™ (FP)	Preincubation Study	Km	1mM	Assay Service	CDK Panel	K192 Panel
GSK3α(GSK3A)	○		○					○	○	○	○	○		
GSK3β(GSK3B)	○	○	○		○			○	○	○	○	○		
Haspin(GSG2)	○		○					○	○	○				
HCK	○		○			○		○		○	○	○		
HER2(ERBB2)	○		○					○		○	○			
HER4(ERBB4)	○		○					○	○	○	○			
HGK(MAP4K4)	○		○	○				○		○	○			
HH498(TNNI3K)	○											○		○
HIPK1	○		○					○	○	○				
HIPK2	○		○					○	○	○		○		○
HIPK3	○		○					○	○	○	○	○		○
HIPK4	○		○					○		○	○	○		○
HPK1(MAP4K1)	○		○					○		○		○		○
HRI(EIF2AK1)	○													
ICK	○											○		○
IGF1R	○	○	○			○		○	○	○	○	○		○
IGF1R [non-activated]		○												
IKKα(CHUK)	○	○		○				■		○				
IKKα(CHUK) [inactive mutant]		○												
IKKβ(IKBKB)	○	○	○					○	○	○	○			
IKKε(IKBKE)	○		○	○				○		○		○		○
INSR	○	○	○					○	○	○	○	○		○
INSR [non-activated]		○												
IRAK1	○			○				■		○		○		
IRAK3												○		○
IRAK4	○	○	○	○				○		○	○	○		○
IRR(INSRR)	○		○					○		○	○			
ITK	○	○	○					○		○	○	○		○
JAK1												○		
JAK1(JH1)	○	○	○			○		○		○	○	○		
JAK2												○		
JAK2 [V617F]												○		○
JAK2(JH1)	○	○	○			○		○	○	○	○	○		
JAK2(JH1JH2)	○		○											
JAK2(JH1JH2) [V617F]	○		○											
JAK3												○		○
JAK3(JH1)	○		○			○		○		○	○			
JNK1(MAPK8)	○							○	○	○	○	○		○
JNK1(MAPK8) [inactive mutant]	○													
JNK1(MAPK8) [inactive]	○													
JNK2(MAPK9)	○							○	○	○	○	○		○
JNK2(MAPK9) [inactive]	○													
JNK3(MAPK10)	○							○	○	○	○	○		○
KDR(VEGFR2)	○	○	○					○	○	○	○			
KHS1(MAP4K5)												○		○
KIT	○	○	○					○		○	○	○		
KIT [A829P]												○		
KIT [D816E]	○		○					○		○	○			
KIT [D816H]												○		
KIT [D816V]	○		○					○		○	○	○		
KIT [D816Y]	○		○					○		○	○			
KIT [L576P]												○		
KIT [T670I/D816V]	○													
KIT [T670I]	○		○					○		○	○			
KIT [V559D,T670I]												○		
KIT [V559D,V654A]												○		
KIT [V559D]												○		
KIT [V560G/D816V]	○		○											
KIT [V560G]	○		○					○		○	○			
KIT [V654A/D816V]	○													
KIT [V654A]			○					○		○	○			
KIT [non-activated]		○												
LATS1												○		○
LATS2			○					○		○	○	○		○
LATS1/MOBKL1A	○													
LATS2/MOBKL1A	○	○												
LCK	○	○	○			○		○	○	○	○	○		○
LIMK1	○											○		○
LIMK2	○	○										○		○
LKB1(STK11)												○		○
LKB1(STK11)/MO25α/STRADα	○													
LOK(STK10)	○		○					○		○		○		○
LRRK2	○											○		○
LRRK2 [G2019S]	○											○		
LRRK2 [I2020T]												○		
LRRK2 [R1441C]												○		
LTK	○		○					○		○	○	○		○

Products & Services by Kinase

<div>Products & Services</div> <div>Kinases name</div>	Products							Services						
	Kinase Proteins	Biotiny-lated Kinases	Kinase Protein Assay Kit					Cell-Free Assays				Cell-Based Assays		
								Mobility Shift Assay/IMAP™		ADP-Glo™ (ATP=Km)	NanoBRET™ TE Intracellular Kinase Cell-Based Assay Services	Assay Service	CDK Panel	K192 Panel
								MSA/IMAP™	ATP conc.					
			MSA	FP (IMAP™)	ELISA	TR-FRET	ADP-Glo™	■: IMAP™ (FP)	Preincubation Study	Km	1mM			
LTK [I565N]	<input type="radio"/>													
LTK [L590M]	<input type="radio"/>													
LTK [L592F]	<input type="radio"/>													
LTK [L650F]	<input type="radio"/>													
LYNa	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>			<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>			
LYNb	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		
LZK(MAP3K13)	<input type="radio"/>				<input type="radio"/>							<input type="radio"/>		
MAP2K1	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>				
MAP2K1 [F129L]	<input type="radio"/>													
MAP2K1 [inactive mutant]	<input type="radio"/>													
MAP2K1 [inactive]	<input type="radio"/>	<input type="radio"/>												
MAP2K1 [P124L]	<input type="radio"/>													
MAP2K2	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>				
MAP2K3	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>				
MAP2K4	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>				
MAP2K5	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		
MAP2K6	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		
MAP2K6 [inactive mutant]	<input type="radio"/>													
MAP2K6 [inactive]	<input type="radio"/>													
MAP2K7	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>				
MAP2K7 [inactive mutant]	<input type="radio"/>													
MAP2K7 [inactive]	<input type="radio"/>													
MAP3K1	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>				
MAP3K2	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MAP3K3	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MAP3K4	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MAP3K5	<input type="radio"/>	<input type="radio"/>						<input type="radio"/>		<input type="radio"/>				
MAP3K6	<input type="radio"/>													
MAP3K14	<input type="radio"/>													
MAP4K2	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MAPKAPK2	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>				<input type="radio"/>	<input type="radio"/>	<input type="radio"/>				
MAPKAPK3	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>				<input type="radio"/>	<input type="radio"/>	<input type="radio"/>				
MAPKAPK5	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>				<input type="radio"/>	<input type="radio"/>	<input type="radio"/>				
MARK1	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>				
MARK2	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MARK3	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>		
MARK4	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>
MAST3												<input type="radio"/>		<input type="radio"/>
MAST4												<input type="radio"/>		<input type="radio"/>
MELK	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>				<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MELK [T460M]												<input type="radio"/>		
MER(MERTK)	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>			<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MER(MERTK) [A708S]												<input type="radio"/>		
MET	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>			<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MET [D1228H]	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>		
MET [D1228N]	<input type="radio"/>	<input type="radio"/>										<input type="radio"/>		
MET [F1200I]												<input type="radio"/>		
MET [M1250T]	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>		
MET [P991S]												<input type="radio"/>		
MET [T1173I]												<input type="radio"/>		
MET [T992I]												<input type="radio"/>		
MET [V1092I]												<input type="radio"/>		
MET [Y1230A]												<input type="radio"/>		
MET [Y1230C]		<input type="radio"/>										<input type="radio"/>		
MET [Y1230D]												<input type="radio"/>		
MET [Y1230H]	<input type="radio"/>	<input type="radio"/>										<input type="radio"/>		
MET [Y1235D]	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>		
MINK(MINK1)	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>		<input type="radio"/>				
MLK1(MAP3K9)	<input type="radio"/>	<input type="radio"/>			<input type="radio"/>			<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MLK2(MAP3K10)	<input type="radio"/>	<input type="radio"/>			<input type="radio"/>			<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MLK3(MAP3K11)	<input type="radio"/>	<input type="radio"/>			<input type="radio"/>			<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		<input type="radio"/>
MLK4	<input type="radio"/>											<input type="radio"/>		<input type="radio"/>
MNK1(MKNK1)	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>				<input type="radio"/>	<input type="radio"/>	<input type="radio"/>				
MNK2(MKNK2)	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>
MOK												<input type="radio"/>		
MOS	<input type="radio"/>				<input type="radio"/>			<input type="radio"/>		<input type="radio"/>				
MRCKα(CDC42BPA)	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>				<input type="radio"/>		<input type="radio"/>				
MRCKβ(CDC42BPB)	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>		<input type="radio"/>				
MSK1(RPS6KA5)	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>				<input type="radio"/>	<input type="radio"/>	<input type="radio"/>				
MSK2(RPS6KA4)	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>				<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>
MSSK1(STK23)	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>		<input type="radio"/>				
MST1(STK4)	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>
MST2(STK3)	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>					<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>
MST3(STK24)	<input type="radio"/>		<input type="radio"/>					<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		
MST4	<input type="radio"/>		<input type="radio"/>	<input type="radio"/>				<input type="radio"/>		<input type="radio"/>		<input type="radio"/>		
MTOR(FRAP)/MLST8	<input type="radio"/>											<input type="radio"/>		
MUSK	<input type="radio"/>	<input type="radio"/>	<input type="radio"/>					<input type="radio"/>		<input type="radio"/>	<input type="radio"/>	<input type="radio"/>		<input type="radio"/>

Products & Services by Kinase

	Products							Services								
								Cell-Free Assays				Cell-Based Assays				
<div>Products & Services</div>	Kinase Proteins	Biotiny-lated Kinases	Kinase Protein Assay Kit					Mobility Shift Assay/ IMAP™		ADP-Glo™ (ATP=Km)	NanoBRET™ TE Intracellular Kinase Cell-Based Assay Services					
								MSA/IMAP™						ATP conc.		
			Kinases name			MSA	FP (IMAP™)	ELISA	TR-FRET		ADP-Glo™	■: IMAP™ (FP)	Preincubation Study	Km	1mM	
MUSK [non-activated]		○														
MYLK3														○		○
MYLK4	○	○												○		○
MYT1(PKMYT1)	○													○		○
NDR1(STK38)	○		○					○		○				○		○
NDR2(STK38L)	○		○					○		○				○		○
NEK1	○		○					○		○	○			○		○
NEK2	○		○					○	○	○	○			○		○
NEK3	○													○		○
NEK4	○		○					○	○	○				○		○
NEK5														○		○
NEK6	○		○					○		○	○			○		
NEK7	○		○					○		○	○					
NEK9	○	○	○					○		○	○			○		○
NEK11	○													○		○
NIM1K(MGC42105)	○		○	○				○	○	○				○		○
NLK	○													○		○
NPM1-ALK	○		○					○	○	○	○					
NRK(NESK)														○		
NuaK1(ARK5)	○		○	○				○	○	○	○			○		○
NuaK2	○	○	○					○	○	○	○			○		
OSR1(OXSR1)	○															
p38α(MAPK14)	○	○	○	○				○	○	○	○			○		○
p38α(MAPK14) [inactive mutant]	○															
p38α(MAPK14) [inactive]	○	○														
p38α(MAPK14) [T106M]														○		
p38β(MAPK11)	○		○	○				○	○	○	○			○		○
p38γ(MAPK12)	○		○	○				○	○	○	○					
p38δ(MAPK13)	○		○	○				○	○	○	○					
p70S6K(RPS6KB1)	○		○	○				○	○	○	○					
p70S6Kβ(RPS6KB2)	○		○	○				○	○	○						
PAK1	○		○					○	○	○	○					
PAK2	○		○					○	○	○	○					
PAK3	○		○													
PAK4	○	○	○					○		○				○		○
PAK5(PAK7)	○		○					○	○	○	○			○		
PAK6	○	○	○					○		○				○		○
PASK	○		○	○				○		○	○					
PBK	○		○	○				○		○	○					
PCTAIRE1(CDK16)/CycY	○													○	○	○
PDGFRα(PDGFRα)	○		○					○	○	○	○					
PDGFRα(PDGFRα) [non-activated]		○														
PDGFRα(PDGFRα) [D842V]	○		○					○	○	○	○					
PDGFRα(PDGFRα) [T674I]	○		○					○		○	○					
PDGFRα(PDGFRα) [V561D]	○		○					○	○	○	○			○		
PDGFRβ(PDGFRβ)	○	○	○			○		○	○	○	○					
PDGFRβ(PDGFRβ) [non-activated]		○														
PDHK1(PDK1)	○															
PDHK2(PDK2)	○		○					○		○						
PDHK3(PDK3)	○															
PDHK4(PDK4)	○		○					○		○						
PDK1(PDPK1)	○							○		○	○					
PEK(EIF2AK3)	○			○				■		○						
PGK(PRKG1)	○		○	○				○		○						
PHKG1	○		○	○				○		○				○		○
PHKG2	○		○					○		○				○		○
PI4KA														○		
PI4KB	○													○		
PIK3C3	○													○		
PIK3CA/PIK3R1	○	○					○					○		○		
PIK3CA [C420R]/PIK3R1														○		
PIK3CA [E542K]/PIK3R1		○					○							○		
PIK3CA [E545A]/PIK3R1														○		
PIK3CA [E545K]/PIK3R1		○					○							○		
PIK3CA [H1047L]/PIK3R1														○		
PIK3CA [H1047R]/PIK3R1		○					○							○		
PIK3CA [H1047Y]/PIK3R1														○		
PIK3CA [I800L]/PIK3R1														○		
PIK3CA [M1043I]/PIK3R1														○		
PIK3CA [P539R]/PIK3R1		○					○							○		
PIK3CA [Q546K]/PIK3R1														○		
PIK3CA [R88Q]/PIK3R1		○					○							○		
PIK3CB/PIK3R1	○	○					○							○		
PIK3CD/PIK3R1	○	○					○							○		
PIK3CG		○														
PIKFYVE(PIP5K3)	○	○					○							○		
PIM1	○		○	○				○	○	○	○					

		Products						Services							
								Cell-Free Assays				Cell-Based Assays			
<div>Products & Services</div>	Kinase Proteins	Biotiny-lated Kinases	Kinase Protein Assay Kit					Mobility Shift Assay/IMAP™			ADP-Glo™ (ATP=Km)	NanoBRET™ TE Intracellular Kinase Cell-Based Assay Services			
								MSA/IMAP™		ATP conc.					
			Kinases name	■: IMAP™ (FP)	Preincubation Study	Km	1mM	Assay Service	CDK Panel	K192 Panel					
PIM2	○		○	○				○		○	○				
PIM3	○		○					○		○	○		○		
PIP4K2A	○						○					○			
PIP4K2B	○						○					○			
PIP4K2C	○												○		
PIP5K1A	○						○					○			
PIP5K1B	○						○					○	○		
PIP5K1C	○						○					○			
PIP5KL1	○						○					○			
PKACα(PRKACA)	○	○	○	○				○		○	○		○		○
PKACβ(PRKACB)	○		○					○		○	○		○		○
PKACγ(PRKACG)	○		○					○		○					
PKCα(PRKCA)	○	○	○	○				○		○	○		○		
PKCβ1(PRKCB1)	○		○	○				○		○			○		
PKCβ2(PRKCB2)	○		○					○		○					
PKCγ(PRKCG)	○	○	○					○		○	○		○		
PKCδ(PRKCD)	○		○	○				○					○		
PKCε(PRKCE)	○		○	○				○	○	○	○		○		○
PKCζ(PRK CZ)	○		○	○				○		○					
PKCη(PRKCH)	○	○	○	○				○		○			○		
PKCθ(PRK CQ)	○		○	○				○		○			○		
PKCι(PRK CI)	○		○	○				○		○					
PKD1(PRK D1)	○		○	○				○	○	○					
PKD2(PRK D2)	○		○	○				○	○	○	○				
PKD3(PRK D3)	○		○	○				○		○					
PKN1	○	○						■		○					
PKN2	○														
PKN3	○														
PKR(EIF2AK2)	○			○				■		○					
PLK1	○	○	○	○				○		○	○		○		
PLK2	○			○				■		○			○		○
PLK3	○		○	○				○		○	○		○		○
PLK4													○		○
PRKX	○		○	○				○		○			○		○
PYK2(PTK2B)	○	○	○					○	○	○	○		○		○
QIK(SNF1LK2)	○		○					○	○	○	○		○		○
RAF1(CRAF)	○	○			○			○		○			○		
RET	○	○	○					○	○	○	○		○		○
RET [G691S]	○	○	○					○	○	○	○				
RET [G691S] [non-activated]		○													
RET [G810C]	○														
RET [G810C/M918T]	○														
RET [G810R]	○														
RET [G810S]	○														
RET [M918T]	○	○	○					○	○	○	○		○		

		Products						Services								
								Cell-Free Assays				Cell-Based Assays				
Products & Services	Kinase Proteins	Biotiny- lated Kinases	Kinase Protein Assay Kit					Mobility Shift Assay/IMAP™			ADP-Glo™ (ATP=Km)	NanoBRET™ TE Intracellular Kinase Cell-Based Assay Services				
								MSA/IMAP™		ATP conc.						
			Kinases name			MSA	FP (IMAP™)	ELISA	TR-FRET	ADP-Glo™		■: IMAP™ (FP)	Preincubation Study	Km	1mM	Assay Service
skMLCK(MYLK2)	○		○						○	○	○	○		○		○
SLK	○		○						○		○			○		○
Smad1	○															
Smad3	○															
smMLCK(MYLK)	○															
SNRK														○		○
SPHK1	○	○	○						○	○	○					
SPHK2	○	○	○						○		○					
SRC	○	○	○						○	○	○	○		○		
SRM(SRMS)	○		○						○	○	○	○		○		○
SRPK1	○	○		○					■		○					
SRPK2	○		○	○					○		○					
SSTK(TSSK6)	○															
STK16														○		○
STK32A														○		
STK32B														○		○
STK33	○													○		○
STK35 (CLIK1)														○		○
STK36														○		○
STLK3(STK39)	○															
SYK	○	○	○			○			○		○	○				
SYK [non-activated]		○														○
TAK1-TAB1(MAP3K7)	○	○	○		○				○		○	○				
TAOK2	○		○						○		○	○				
TAOK3	○															
TBK1	○	○	○						○	○	○	○		○		○
TEC	○		○						○	○	○	○		○		○
TESK1	○													○		
TGFβR1(TGFBR1)(ALK5)	○	○				○								○		
TGFβR2(TGFBR2)	○													○		
TIE1														○		○
TIE2(TEK)	○	○	○			○			○	○	○	○		○		○
TIE2(TEK) [A1124V]														○		
TIE2(TEK) [P883A]														○		
TIE2(TEK) [R849W]														○		
TIE2(TEK) [Y1108F]														○		
TIE2(TEK) [Y897C]														○		
TIE2(TEK) [Y897S]														○		
TLK1	○													○		○
TLK2	○													○		○
TNIK	○	○	○	○					○		○	○				
TNK1	○		○						○		○	○		○		○
TRKA(NTRK1)	○	○	○			○			○		○	○		○		○
TRKA(NTRK1) [non-activated]		○														

Products & Services by Kinase

		Products						Services						
								Cell-Free Assays				Cell-Based Assays		
Products & Services <														



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

Carna's wholly-owned subsidiary

CarnaBio USA, Inc.

209 West Central Street, Suite 127, Natick, MA 01760 USA

TEL: +1-508-650-1244

Toll-Free: +1-888-645-1233

E-mail: orders@carnabio.com / FAX: +1-508-650-1722