



# KINASE RESEARCH TOOLS

FOR CELL SIGNALING DISCOVERY

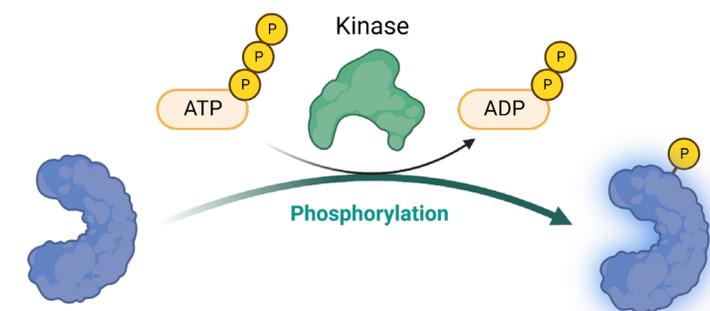
Active Proteins | Assay Kits | Substrates | Inhibitors  
CRISPR/Cas9 Kinase Knock-out Lentivirus Library



# A Master Controller of Cell Functions

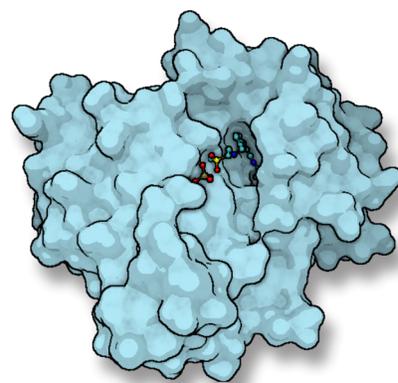
Protein phosphorylation regulates a large number of cellular processes including cellular division, metabolism, signal transduction, and many more. There are over 500 kinases encoded by the human genome and it has been estimated that kinases regulate approximately 50% of cellular functions. Kinase is the name given to an enzyme that catalyzes the transfer of a phosphate group from ATP to another molecule. Kinases can be broadly separated by their substrate into protein, lipid, nucleotide, and carbohydrate kinases with the protein kinases being further classified based on their cellular location and the specific amino acid(s) they phosphorylate. They are now key therapeutic targets across many disease areas.

## Phosphorylation



Target protein

Images created with Biorender.com.



EGFR kinase domain complexed with an inhibitor, Lapatinib. (PDB ID: 1XKK)

## Our Advantages



### Produced In-House

- Made in the USA at our San Diego, CA laboratory
- Get customized, personal support directly from the source



### Committed to Excellence

- ISO 9001:2015-certified Quality Management System
- Lot-specific quality testing



### Protein Expression & Screening Services

- Unique panel of 80+ assays for evaluation of lead compounds
- Select from IC<sub>50</sub> determination or single point concentrations
- Perform follow-up studies using the same proteins manufactured in-house



### Expansive Portfolio

- Active kinase complexes: mTOR, 10+ PI3 Kinase complexes, AMPK, and more
- Active kinase mutants: TRKC, LRRK2, JAK2, cKit, and 35+ EGFR mutants
- 100+ activity assay kits
- CRISPR/Cas9 Kinase Knockout Lentiviruses

# Bioactive Recombinant Proteins

BPS Bioscience specializes in manufacturing bioactive enzymes and protein complexes, offering over 2000 protein products critical for drug discovery. Our kinase portfolio consists of proteins from six kinase families: lipid, receptor tyrosine, serine-threonine, cytoplasmic tyrosine, pyruvate, and fructokinase. Our offerings include wild-type and mutant variants of key proteins, including JAK2, TRKC, LRRK2, FGFR3, and EGFR.

Protein substrates and inhibitors are also available.

Protein options include:



## HiP™ Proteins (High Purity, Low Aggregation)

- >90% purity
- Verified via HPLC to guarantee <10% aggregation
- Ideal for studying enzyme activity, ligand binding, protein folding, and more

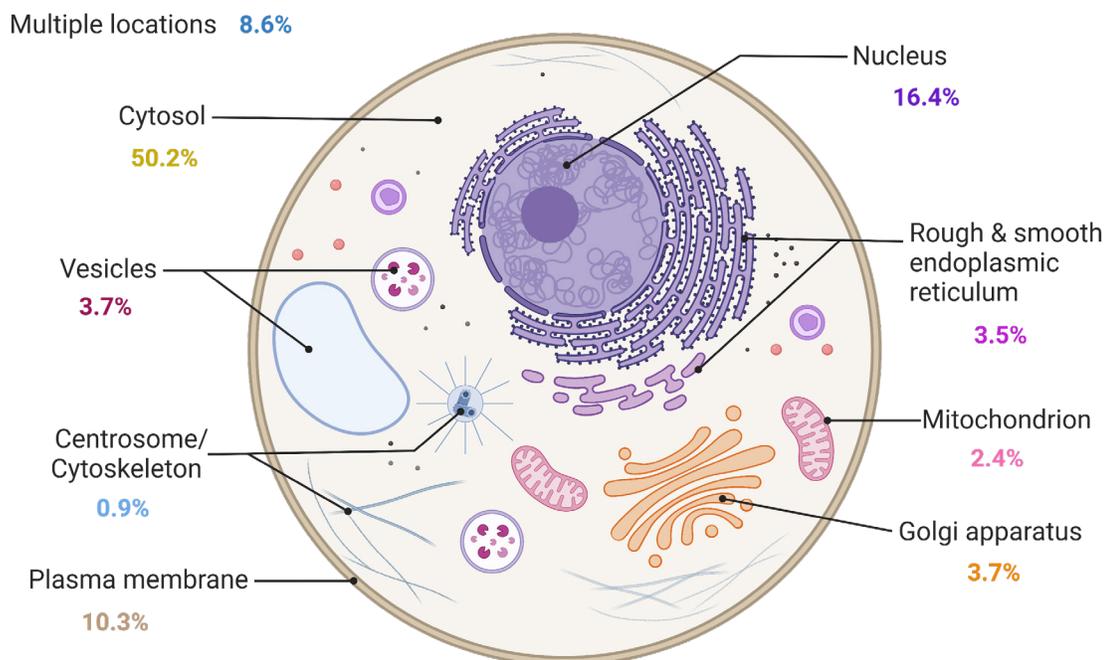
## Bulk Production and Customization

Product sizing, formulations, tags, mutations, and activity testing can be customized to meet your research needs.

## Featuring

- Active complexes, including AMPK, PI3 Kinase, CDK/Cyclin, and more
- Active protein domains, including JAK2 JH2, JAK3 JH2, TYK2 JH2, PLK1 Polo Box, and more.

## Where are kinases found in the cell?



Data from: Zhang, H. et al. 2021. eLife. [64943](https://doi.org/10.7554/eLife.64943). Image created with Biorender.com.

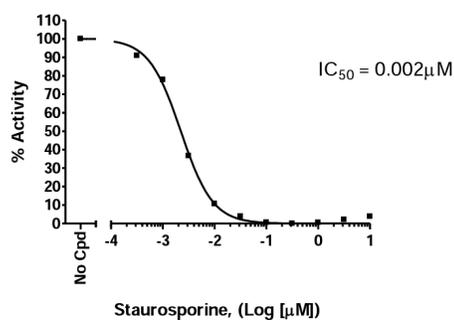
# Biochemical Assay Kits

Our 100+ kinase assay kits enable your compound and biotherapeutic drug discovery by providing specific, highly sensitive analysis in a variety of formats. Many of our kits are provided as homogenous assays, with no wash steps, for efficient high throughput screening.

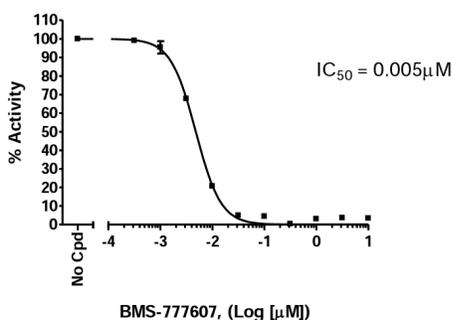
Compare assay formats:

AlphaLISA™	Chemiluminescent	Colorimetric	Fluorescence Polarization	Fluorogenic	Luminescent	TR-FRET	UV Spectrophotometric		
✓	✓	✓	✓	✓	✓	✓	✓		High specificity
✓	✓		✓	✓	✓	✓	✓		High sensitivity with low background noise.
✓			✓	✓	✓	✓	✓		Simple operation
✓			✓	✓	✓	✓	✓		Homogeneous, no-wash format makes it simple for high throughput screening.
✓			✓	✓		✓	✓		Real-time readings do not require equilibrium conditions.
✓			✓			✓	✓		Small volume reactions save on sample consumption.
✓	✓			✓					Large dynamic range, up to 4-5 logs.
			✓			✓			Highly stable signal allows for repetitive measurements.
		✓					✓		Common, low-cost instrumentation
						✓			Use of lanthanide donor fluorophores minimizes photobleaching.

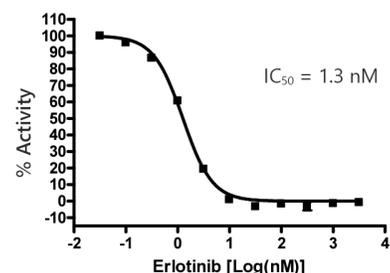
JAK2 Inhibition by Staurosporine



Tyro3 Inhibition by BMS-777607



EGFR Inhibition by Erlotinib



# Biochemical Screening Services

BPS Bioscience offers rapid and dependable kinase screening services to determine  $IC_{50}$  values of small molecules, peptides, and antibodies against requested targets. Our team of experts along with our broad services portfolio make it easy to:

- Screen for inhibitors/targets
- Select from  $IC_{50}$  determination and single point concentrations
- Perform follow-up studies using the same proteins manufactured in-house
- Get questions answered or project guidance in a time-efficient manner

## Our Deliverables to You

### Detailed Results



Extensive report with raw and analyzed data, graphs, and detailed protocols. Typically includes positive controls for inhibition.

### Fast Turnaround



Receive results within weeks after our receipt of your compounds.

### Quality & Consistency



Proteins and enzymes synthesized and QC'd in-house to ensure the highest level of inter- and intra-assay consistency.

### Extensive Selection



Choose from over 80 in-house developed kinase assays, including many novel assay kits.

### Platform Options



Choose the format that works best for your experiment or perform orthogonal testing on multiple platforms to verify your results.

## Why choose BPS Bioscience for your screening projects?

- We can provide end-to-end customization from protein production to assay development, completing with compound screening.
- We have extensive expertise and experience in developing and manufacturing custom and off-the-shelf kinase proteins and assay kit products.
- We have helped accelerate projects across large pharma, vaccine developers, biotech and basic research institutions.
- Our high quality projects have returned excellent customer satisfaction scores.

*Give us a try today.*

# CRISPR/Cas9 Kinase Knockout Library

Our CRISPR/Cas9 Kinase Knockout Lentivirus Library consists of ready-to-use, VSV-G pseudotyped integrating lentiviruses intended for use in generating kinase knockouts for screening purposes.

## Array Format (#78487)



Targets 619 human kinases and pseudokinases



5 sgRNAs per gene



150 control sgRNAs included



Puromycin selection

- 649 vials come with 200  $\mu$ l of lentivirus each at a titer  $\geq 1 \times 10^7$  TU/ml; the exact titer is provided with each shipment.

## Advantages

- All-in-One: each vial contains lentiviral particles to transduce the Cas9 gene, 5 sgRNAs targeting the kinase of interest, and a puromycin selection marker. There is no requirement for the target cell to already express Cas9.
- Safe: requires only Biosafety Level 2 (BSL-2).
- Each sgRNA lentiviral particle is individually constructed, sequence-verified, individually cultured, and titered to ensure high quality and representation across the entire library.
- VSV-G pseudoviruses transduce most mammalian cells, including primary and non-dividing cells.

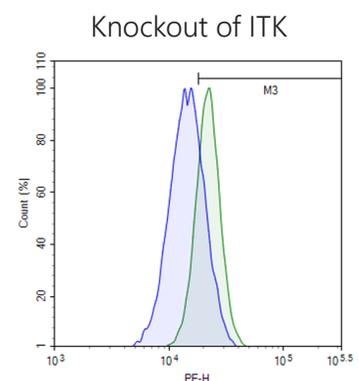
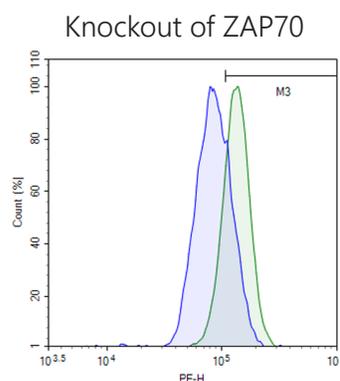
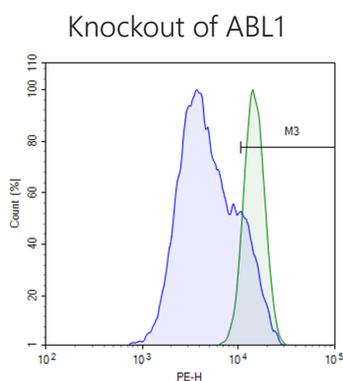
## Custom Panels

- The Custom Panel option allows you to pick between 2-50 kinases out of the full human kinase library. One vial of 200  $\mu$ l of lentivirus is provided for each target.
- Download the full target list and request a quote here: [bpsbioscience.com/crispr-kinasepanel](https://bpsbioscience.com/crispr-kinasepanel)

## Single Target (#78488)

- If you are only interested in a single kinase target to knock out, this option is available.
- Receive two vials, each with 200  $\mu$ l of lentivirus at a titer  $\geq 1 \times 10^7$  TU/ml.

Flow cytometry confirmation of kinase knockouts in Jurkat cells:



Assay Kit Components	Catalog#	Biochemical Assay Kits	Catalog#
ATP (500 μM)	79686	Chemi-Verse™ CDK3/CyclinE1 Kinase Assay Kit	78884
		Chemi-Verse™ CDK5/p35 Kinase Assay Kit	82249
		Chemi-Verse™ CDK8/Cyclin C Kinase Assay Kit	78886
		Chemi-Verse™ CDK9/CyclinK Kinase Assay Kit	82237
		Chemi-Verse™ CDK9/CyclinT Kinase Assay Kit	82233
		Chemi-Verse™ CLK1 Kinase Assay Kit	82145
		Chemi-Verse™ CLK2 Kinase Assay Kit	82146
		Chemi-Verse™ CLK3 Kinase Assay Kit	82150
		Chemi-Verse™ ERK2 Kinase Assay Kit	78899
		Chemi-Verse™ Kinase Aurora C Assay Kit	82096
		Chemi-Verse™ LYN A Kinase Assay Kit	82224
		Chemi-Verse™ MAPKAPK2 Kinase Assay Kit	82129
		Chemi-Verse™ MAPKAPK5 Kinase Assay Kit	82236
		Chemi-Verse™ MEK1 Kinase Assay Kit	82100
		Chemi-Verse™ NUA1 Kinase Assay Kit	82231
		Chemi-Verse™ NUA2 Kinase Assay Kit	82232
		Chemi-Verse™ PAK2 Kinase Assay Kit	82221
		Chemi-Verse™ PAK3 Kinase Assay Kit	82222
		Chemi-Verse™ PAK4 Kinase Assay Kit	82181
		Chemi-Verse™ PAK6 Kinase Assay Kit	82182
		Chemi-Verse™ PAK7 Kinase Assay Kit	82183
		Chemi-Verse™ PI3 Kinase (p110α/p85α) Kinase Assay Kit	82108
		Chemi-Verse™ PI3 Kinase (p120γ) Kinase Assay Kit	82104
		Chemi-Verse™ PI3 Kinase p110α (E542K)/p85α Kinase Assay Kit	82113
		Chemi-Verse™ PI3 Kinase p110α (E545K)/p85α Kinase Assay Kit	82109
		Chemi-Verse™ PI3 Kinase P110α (H1047R)/P85α Kinase Assay Kit	82103
		Chemi-Verse™ PI3 Kinase P110α(H1047L)/P85α Kinase Assay Kit	82120
		Chemi-Verse™ SIK1 Kinase Assay Kit	82105
		Chemi-Verse™ SIK2 Kinase Assay Kit	82106
		Chemi-Verse™ TNK1 Kinase Assay Kit	82159
		Chemi-Verse™ TTBK1 Kinase Assay Kit	82162
		Chemi-Verse™ TTBK2 Kinase Assay Kit	82225
		Chemi-Verse™ VRK1 Kinase Assay Kit	82230
		CHK1 Kinase Assay Kit	78418
		CHK2 Kinase Assay Kit	78421
		cKIT (D816V) Assay Kit	79889
		cKIT (V654A) Assay Kit	79890
		cKIT Assay Kit	79815

Biochemical Assay Kits	Catalog#	Biochemical Assay Kits	Catalog#
CSF1R Kinase Assay Kit	78829	JAK2-JH2 Pseudokinase Domain (V617F, W659A, W777A, F794H) Inhibitor Screening Assay Kit	78209
CSK Kinase Assay Kit	78818	JAK2-JH2 Pseudokinase Domain (W659A, W777A, F794H) Inhibitor Screening Assay Kit	78208
DYRK1B Kinase Assay Kit	78401	JAK3 (Janus Kinase 3) Assay Kit	79521
DYRK2 Kinase Assay Kit	78392	LCK Assay Kit	79794
EGFR (T790M, C797S) (del746-750) Kinase Assay Kit	78595	Lumi-Verse™ VEGFR2 (KDR) Kinase Assay Kit	78857
EGFR (T790M/C797S/ L858R) Kinase Assay Kit	40326	MAP3K14 Kinase Assay Kit	78440
EGFR Kinase Assay Kit	40321	MNK1 Kinase Assay Kit	78032
EGFR(L858R) Kinase Assay Kit	40324	NEK6 Kinase Assay Kit	79992
EGFR(T790M) Kinase Assay Kit	40323	NEK7 Kinase Assay Kit	78850
EGFR(T790M/L858R) Kinase Assay Kit	40322	p70S6K (S6K1) Kinase Assay Kit	78806
EIF2AK2 Kinase Assay Kit	78433	PDGFRα (D842I) Assay Kit	79761
EPHA2 Assay Kit	78002	PDGFRα (D842V) Assay Kit	79828
EPHA4 Kinase Assay Kit	78592	PDGFRα (D842Y) Assay Kit	79760
EPHB4 Kinase Assay Kit	78009	PDK3 Kinase Assay Kit	78286
FAK Kinase Assay Kit	40722	PI3Kα (p110α/p85) Assay Kit	79781
FGFR2 Assay Kit	79804	PI3Kβ (p110β/p85α) Assay Kit	79802
FGFR3 (L496V) Assay Kit	79818	PI3Kγ (p110γ/PIK3R5) Assay Kit	79803
FGFR3 (V443L) Assay Kit	79819	PI3Kδ (p110δ/p85α) Assay Kit	79799
FGFR3 (V443M) Assay Kit	79817	PIM1 Assay Kit	79885
FGFR4 Kinase Assay Kit	78816	PIM2 Assay Kit	79883
FGR Kinase Assay Kit	82292	PIM3 Assay Kit	79886
FLT1 Kinase Assay Kit	78019	Rat ITK Kinase Assay Kit	78513
FLT3 Kinase Assay Kit	79797	RET Assay Kit	79566
FYN Assay Kit	78003	RIPK1 Kinase (Mouse) Assay Kit	79897
GSK3α Assay Kit	79698	RIPK1 Kinase Assay Kit	79560
GSK3β Assay Kit	79700	RIPK1 Kinase Assay Kit (384-well)	79689
HER2 Kinase Assay Kit	40721	RIPK2 Kinase Assay Kit	79737
HER4 Assay Kit	78005	ROCK1 Kinase Assay Kit	78386
HIPK3 Kinase Assay Kit	78382	ROCK2 Kinase Assay Kit	78407
HPK1 Assay Kit	79775	ROS1 Kinase Assay Kit	78189
IKKβ Kinase Assay Kit	78507	SPHK1 Assay Kit	78026
IRAK1 Kinase Assay Kit	78523	SRC Assay Kit	79680
IRAK4 Kinase Assay Kit	78524	STK3 (MST2) Kinase Assay Kit	82140
ITK Kinase Assay Kit	78429	STK4(MST1) Kinase Assay Kit	78500
JAK1 (Janus Kinase 1) Assay Kit	79518	SYK Assay Kit	79671
JAK1 (JH2 Pseudokinase Domain) Inhibitor Screening Assay Kit	78891	TAK1-TAB1 Kinase Assay Kit	78037
JAK2 (Janus Kinase 2) Assay Kit	79520	TBK1 Assay Kit	79871
JAK2 JH2 Pseudokinase Domain Inhibitor Screening Assay Kit	78111	TGFβR1 (ALK5) Kinase Assay Kit	78819

Biochemical Assay Kits		Catalog#	Inhibitors/Activators		Catalog#
Tie2 Kinase Assay Kit		78573	CHIR-99021 (CT99021) HCl		27615-3
TNIK Kinase Assay Kit		78370	Doramapimod		27034
TrkA Assay Kit		79548	Fasudil		27029
TRKB Kinase Assay Kit		78591	Forskolin		27067
TRKC Kinase Assay Kit		78528	Forskolin		27068
TTK (MPS-1) Kinase Assay Kit		78356	GDC-0941		27000
TYK2 (Tyrosine Kinase 2) Assay Kit		79527	Gefitinib		27032
TYK2 JH2 Pseudokinase Domain Inhibitor Screening Assay Kit		78107	Geldanamycin		27008
TYR Phosphorylation Assay Kit		79502	HA-1077		27030
TYRO3 Kinase Assay Kit		79593	IBT6A		82075
ULK1 Kinase Assay Kit		78362	K252a		27033
ULK2 Kinase Assay Kit		78367	I-[(3R)-3-[4-amino-3-(4-phenoxyphenyl)-1Hpyrazolo [3,4-d]pyrimidin-1-yl]-1-piperidinyl]-2-propyn-1-one		82071
VEGFR2 (KDR) Kinase Assay Kit		40325	LY-294002		27038
VEGFR3(FLT4) Kinase Assay Kit		79738	MK-1775		82196
Wee1 Kinase Assay Kit		79909	PD 173074		27056
YES1 Assay Kit		79681	PI-103		27502-1
ZAP70 Kinase Assay Kit		78427	PI-103		27502-2
			PP242		27063
			Rapamycin		27062
			Roscovitine		27040
			Ruxolitinib Phosphate		27059
			SGI-1776 free base		27651-1
			SGI-1776 free base		27651-2
			SGI-1776 free base		27651-3
			SMI-4a		27774-1
			SMI-4a		27774-2
			Sorafenib Tosylate		27014
			Staurosporine		27002
			Temsirolimus		27024
			Tofacitinib		27037
			U0126		27012
			Wortmannin		27031
Buffers		Catalog#	Lentiviruses		Catalog#
Kinase Buffer 1		79334	CRISPR/Cas9 Kinase Knockout Lentivirus Library (Array Format)		78487
			Kinase (Human) CRISPR/Cas9 Lentivirus (Integrating)		78488
Cell Lines		Catalog#			
CSF1R / SRE - Reporter HEK293 Recombinant Cell Line		79380			
Human EGFRvIII CHO K1 Cell Line (High or Low Expression)		78145			
TrkA / SRE Reporter - HEK293 Recombinant Cell Line		79798			
VEGFR2/NFAT Reporter HEK293 Recombinant Cell Line		79387			
Inhibitors/Activators		Catalog#			
4-(4-(4-(3-fluorobenzyloxy)-3-chlorophenylamino)quinazolin-6-yl)phenol		82072			
A-769662		27057			
Afatinib		27009			
Axitinib		27064			
BAY-11-7082		27229			
BI-2536		82062			
Cediranib		27036			
CHIR-99021 (CT99021)		27614-2			
CHIR-99021 (CT99021)		27614-3			
CHIR-99021 (CT99021) HCl		27615-2			

Proteins	Catalog#	Proteins	Catalog#
14-3-3 alpha/beta Protein, His-tag (Human)	79023	ALK2 (ACVR1) (Q207D), GST-Tag	101936
14-3-3 alpha/beta Protein, His-tag (Mouse)	79024	ALK2 (ACVR1) (Q207E), GST-Tag	101937
14-3-3 epsilon Protein, His-tag (Human)	79025	ALK2 (ACVR1) (R202I), GST-Tag	101938
14-3-3 epsilon Protein, His-tag (Mouse)	79026	ALK2 (ACVR1) (R206H), GST-Tag (Mouse)	101940
14-3-3 sigma Protein, His-tag (Human)	79037	ALK2 (ACVR1) (R206H), GST-Tag	100714
14-3-3 theta Protein, His-tag (Human)	79038	ALK2 (ACVR1) (R258G), GST-Tag	101941
14-3-3 zeta Protein, His-tag (Human)	79039	ALK2 (ACVR1), FLAG-Tag, Avi-Tag, Biotin-Labeled	100176
14-3-3 zeta Protein, His-tag (Mouse)	79040	ALK2 (ACVR1), GST-tag	40019
Abl Kinase Mutant (T315I), His-tag	40415	ALK3 (BMPR1A) (Q233D), GST-Tag	101942
Abl Kinase, His-tag	40411	ALK3 (BMPR1A) (R486Q), GST-Tag	101943
ABL2, His-tag	40417	ALK4 (ACV1B), GST-tag	40020
ACK1, GST-tag	40400	AMPK (A1/B1/G1), His-tags	40025
AKT1, Active	40003	AMPK (A1/B1/G2), His-tag	40701
AKT1, Inactive, His-tag	40000	AMPK (A1/B1/G2), His-tags	40021
AKT2, Active, His-tag	40011	AMPK (A1/B1/G3), His-tag	40702
AKT2, Inactive, His-tag	40001	AMPK (A1/B1/G3), His-tags	40022
AKT3, His-tag	40012	AMPK (A1/B2/G1), His-tag	40703
ALK (C1156Y), GST-Tag	101915	AMPK (A1/B2/G1), His-tags	40023
ALK (F1174L), GST-Tag	101918	AMPK (A2/ B1/G1), His-tag	40704
ALK (F1174S), GST-Tag	101919	AMPK (A2/B1/G1), His-tags	40024
ALK (G1202R), GST-Tag	101920	AMPK (A2/B2/G1), His-tag	40705
ALK (G1269A), GST-Tag	101921	AMPK (A2/B2/G1), His-tags	40026
ALK (G1269S), GST-Tag	101922	AMPK (A2/B2/G2), His-tag	40706
ALK (I1171N, D1203N), GST-Tag	101923	AMPK (A2/B2/G2), His-tags	40027
ALK (L1152R), GST-Tag	101924	aRAF, GST-Tag, His-Tag	40010
ALK (L1196M), GST-Tag	101926	Ask1, GST-tag	40028
ALK (L1196M, G1202R), GST-Tag	101925	ATF2, GST-Tag	40520
ALK (R1275Q), GST-Tag	101927	ATM, GST-tag	101486
ALK (S1206R), GST-Tag	101928	ATR, GST-Tag	101485
ALK (T1151 L1152insT), GST-Tag	101930	Aurora A Protein Kinase, His-Tag (HEK293-derived)	100112
ALK (T1151M), GST-Tag	101929	Aurora A Protein Kinase, His-Tag (Sf9-derived)	40004
ALK , GST-tag	40214	AURORA C, GST-tag	40178
ALK1 (ACVRL1), GST-tag	40018	Aurora Kinase B, GST-Tag, His-Tag	40002
ALK2 (ACVR1) (G328E), GST-Tag	101931	AXL, FLAG-Tag	100174
ALK2 (ACVR1) (G328V), GST-Tag	101932	Axl, GST-Tag	40180
ALK2 (ACVR1) (G328W), GST-Tag	101933	BCKDK, GST-Tag	101115
ALK2 (ACVR1) (G356D), GST-Tag	101934	BLK, GST-tag	40401
ALK2 (ACVR1) (PF197_198L), GST-Tag	101935	BMX, GST-tag	40402

Proteins	Catalog#	Proteins	Catalog#
BRAF (V600E), GST-His-Tag	40533	CDK12/Cyclin K, GST-Tags	101235
BRAF, GST-tag	40065	CDK13/Cyclin K, GST-Tags	101128
BRAF, His-Tag (High Expression)	41237	CDK14/Cyclin Y, GST-Tag	100602
BRAF/p50, FLAG-tag	40005	CDK16/Cyclin Y, GST-tags	100604
BRK, GST-tag	40403	CDK17/Cyclin Y, GST-tags	100606
BRSK2, His-tag	40016	CDK18/Cyclin Y, GST-tags	100608
BTK(C481F), His-Tag	101040	CDK19/Cyclin C, GST-tags	100593
BTK(C481R), His-Tag	101041	CDK2 (no tag)/CyclinA2, His-GST-tags	41101
BTK(C481S), His-Tag	101039	CDK2/CyclinA2, GST-tag	40101
BTK(P190K), His-Tag	101042	CDK2/CyclinE1, GST-Tag	40102
BTK(T474I), His-Tag	101045	CDK20, GST-Tag (Mouse)	100609
BTK(Y485F), His-Tag	101046	CDK3/CyclinE1, GST-tag	40103
BTK, GST-His-Tag	40405	CDK4(EE,T172A)/Cyclin D1, His-tag	40094
c-Mer, GST-tag (Human)	40254	CDK4, FLAG-Tag	100052
c-MET (del 963-1009), GST-tag	100643	CDK4/CyclinD3, GST-His-Tag	40104
c-Met, GST-tag	40255	CDK5/p25, GST-tag	40105
c-Myc, His-tag	40453	CDK5/p35, GST-tag	40095
c-RAF(Y340D, Y341D), GST-Tag	40008	CDK6, His-Tag	100031
Calmodulin, His-tag	81152	CDK6/CyclinD1, His-tag, GST-tag	40097
CAMK1, Mouse, GST-tag	40030	CDK6/CyclinD3, His-tags	40206
CAMK1 $\beta$ , Mouse, GST-tag	40031	CDK7/CyclinH1/MNAT1, His-tags	40098
CAMK1 $\gamma$ , GST-tag	40032	CDK8/Cyclin C, GST-tags	100433
CAMK1 $\gamma$ , GST-tag	40036	CDK9/CyclinK, GST-tag	40106
CAMK1 $\delta$ , GST-tag	40037	CDK9/CyclinT1, GST-Tag	40307
CAMK2 $\alpha$ , GST-tag	40038	Ceramide Kinase, FLAG-tag	40617
CAMK2 $\beta$ , His-tag	40041	CERK, FLAG-Tag (Mouse)	79413
CAMK2 $\gamma$ , GST-tag	40042	CERK, FLAG-Tag (Rat)	79414
CAMK2 $\delta$ , GST-tag	41109	CHK1, GST-tag	40039
CAMK2 $\delta$ , GST-tag	40043	CHK2, His-tag	40040
CAMK4, GST-tag	40029	CK1 alpha, GST-Tag	79346
CAMKK1, GST-tag	40044	CK1 delta, GST-Tag	79347
CAMKK2, GST-tag	40046	CK1 epsilon, GST-Tag	79349
Casein Kinase 1 $\alpha$ 1	40310	CK1 gamma, GST-Tag	79348
CDC7/DBF4 Complex	100538	CK2a1, GST-tag	40045
CDK1/CyclinA2, GST-Tag	40100	cKit (D816V), His-tag	40252
CDK1/CyclinB1, GST-Tag	40454	cKit (V560G), His-tag	40253
CDK11A/CyclinD3, GST-tags	100595	cKit (V654A), His-tag	40251
CDK12 (696-1082)/CyclinK, GST-Tags,	100998	cKit, His-tag	40250

Proteins	Catalog#	Proteins	Catalog#
CLK1, GST-tag	40196	EGFR (D770GY), GST-Tag	79442
CLK2, GST-tag	40197	EGFR (G719C), GST-Tag	79445
CLK3, GST-tag	40198	EGFR (G719D), GST-Tag	79446
Cofilin 1, GST-tag	40525	EGFR (G719S), GST-Tag	79447
Cot (MAP3K8), GST-His-Tags	40050	EGFR (L747S), GST-Tag	79450
CSF1R, Fc-Fusion, Avi-Tag (Human) HiP™	72550	EGFR (L858R), GST-Tag, His-Tag	40189
CSF1R, Fc-Fusion, Avi-Tag, Biotin-Labeled	79343	EGFR (L861Q), GST-Tag	79451
CSF1R, GST-His-Tag	40227	EGFR (P753S), DEL 747-752, GST-Tag	79439
CSF1R, untagged (FMS, CD115)	79110	EGFR (T790M C797S), DEL 746-750, GST-Tag	79434
CSK, GST-His-Th-Tag	40410	EGFR (T790M C797S), GST-Tag	79453
DAPK1, GST-tag	40051	EGFR (T790M L858R), DEL 746-750, GST-Tag	79435
DAPK2, GST-tag	40049	EGFR (T790M L978I), DEL 746-750, GST-Tag	79436
DAPK3, GST-tag	40052	EGFR (T790M), His-tag, GST-tag	40188
DCAMKL1, GST-tag	40047	EGFR (T790M, C797S, L858R), His-Tag, GST-tag	40351
DCAMKL2, GST-tag	40048	EGFR (T790M, L858R), GST-tag, His-tag	40350
DDR1, Fc Fusion (IgG1), Avi-Tag HiP™	101375	EGFR (V948R), His-Tag	101644
DDR1, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101378	EGFR, DEL 747-749, GST-Tag	79438
DDR1, GST-tag	40184	EGFR, DEL 752-759, GST-Tag	79440
DDR2, Fc Fusion (IgG1), Avi-Tag	101376	EGFR, His-tag, GST-tag	40187
DDR2, Fc Fusion (IgG1), Avi-Tag, Biotin Labeled	101377	EGFR, INS A763-FHEA-Y764, GST-Tag	79427
DDR2, GST-tag	40185	EGFR, INS D770-GF-N771, GST-Tag	79443
DGKa, His-Avi-Tag, Biotin-Labeled	100964	EGFR, INS D770-NPG-N771, GST-Tag	79444
DLK, GST-Tag	79345	EGFR, INS H773-AH-V774, GST-Tag	79448
DMPK, GST-tag	40054	EGFR, INS P772-NPH-H773, GST-Tag	79452
DRAK2	71225	EIF2AK2, GST-tag	40058
DRAK2, GST-tag (Mouse)	100444	EIF2AK3, GST-tag	40059
DRAK2, GST-Tag	71224	eIF4E, GST-tag	40530
DYRK1B, GST-tag	100646	EIF4EBP1, His-tag	40510
DYRK2, His-Tag	40053	EPHA1, GST-tag	40191
EEF2K, GST-tag	40057	EPHA2, His-tag	40190
EGFR (A750P), DEL 747-749, GST-Tag	79437	EPHA3, GST-tag	40192
EGFR (C775S T790M L858R), DEL 746-750, GST-Tag	79431	EPHA4, GST-tag	40193
EGFR (C775S T790M L858R), GST-Tag	79428	EPHA6, GST-tag	40194
EGFR (C797A), DEL 746-750, GST-Tag	79432	EPHB2, His-tag	40200
EGFR (C797A), GST-Tag	79454	EPHB3, GST-tag	40186
EGFR (C797S) GST-Tag	79430	EPHB4, His-tag	40201
EGFR (C797S), DEL 746-750, GST-Tag	79433	EPOR-(GSSG)-6-JAK2(V617F, Y1007F, Y1008F)	100713
EGFR (D761Y), GST-Tag	79441	ERK1	40055

Proteins	Catalog#	Proteins	Catalog#
ERK2, GST-Tag	40299	GRK7, GST-tag	40108
ERK3, GST-Tag	100366	GSK3 $\alpha$	40006
ERK4, GST-Tag	100363	GSK3 $\beta$ , GST-tag, His-tag	40007
ERK7, GST-Tag (Mouse)	100365	GSK $\alpha$ , GST-Tag, His-tag	100814
ERN1 (IRE1), GST-tag	40306	HCK, GST-tag	40440
FAK, GST-Tag	40420	HER2, GST-Tag	40230
FES, His-FLAG-tag	40425	HER3 (E928G), GST-Tag	101645
FGFR1 (V561M), GST-tag	40209	HER3, GST-Tag	101646
FGFR1(FLT2), GST-tag	40210	HER3, His-Tag	101647
FGFR1, Avi-His-Tag, Biotin-labeled (Dephosphorylated)	100692	HER4, GST-tag	40232
FGFR1a (IIIc), Fc Fusion, Avi-Tag	101836	HGK (MAP4K4 ), GST-tag	40109
FGFR2, DEL (429-430), His-Tag (Dephosphorylated)	100356	HIPK1, GST-tag	40179
FGFR2, GST-tag	40211	HIPK3, His-tag	40110
FGFR2C, Fc Fusion, Avi-Tag	101877	HIPK4, GST-tag	40111
FGFR3 (397-806(end)) (K650M) His-Tag (Dephosphorylated)	100909	Histone H3, Full Length, His-tag	52023
FGFR3 (397-806(end)), His-Avi-Tag, Biotin-labeled	100918	Histone H3, Full Length, His-tag	52027
FGFR3 (447-761) (K650M), His-Tag (Dephosphorylated)	100783	HPK1, GST-Tag, His-Avi-Tag (Human)	40398
FGFR3 (447-761), His-Avi-Tag, Biotin-labeled	100613	Human IRE1, His-Avi-tag	40709
FGFR3 (CD333), GST-tag	40212	Human/Mouse ERK2 (K52R), Inactive, His-tag	40056
FGFR3 (K650E), His-Tag (Dephosphorylated)	100782	ICK, GST-tag	40112
FGFR3 (L496V), His-Tag	100142	IGF1R, His-Tag	101217
FGFR3 (N540K) His-Tag (Dephosphorylated)	100907	IKK $\beta$ , GST-Tag	40304
FGFR3 (N540K) His-Tag (Dephosphorylated), Biotin-Labeled	100921	Insulin receptor (INSR), GST-tag	40241
FGFR3 (V443L), His-Tag	100138	IP6K2, Biotin-labeled, His-tag, Avi-tag	40643
FGFR3 (V443M), His-Tag	100140	IP6K2, FLAG-His-tags	40642
FGFR3 (V555F), His-Tag (Dephosphorylated)	100657	IRAK1, His-Tag	40202
FGFR4, GST-tag	40213	IRAK2, GST-tag	40087
FGR, GST-tag	40422	IRAK4, GST-tag	40064
FLT1, His-tag	40223	IRR, GST-tag	40242
Flt3, His-tag	40225	ITK, GST-tag	40445
FRK, GST-tag	40424	JAK1 (JH2 Domain), His-Tag	101946
Fucokinase, FLAG-tag	71077	JAK1, GST-tag	40449
FYN, FLAG-Tag	40433	JAK2 (JH1 domain), His-Tag	40450
Galactokinase, His-tag	50275	JAK2 (JH1, JH2 domain), GST-Tag, Avi-Tag, Biotin (Dephosphorylated)	79495
GCK (MAP4K2), GST-tag	40107	JAK2 (JH1, JH2 domain), His-GST-Tags	40451
GRK2, GST-tag	40060	JAK2 (JH2 Domain) (E596A, V617F, W659A, W777A, F794H), His-Tag	100122
GRK5, GST-tag	40063	JAK2 (JH2 Domain) (V617F, W659A, W777A, F794H), His-Tag	100124
GRK6, GST-tag	40066	JAK2 (JH2 Domain) (W659A, W777A,F794H), His-Tag	100123

Proteins	Catalog#
JAK2 (JH2 Domain), His-Avi-Tag, Biotin-Labeled	79074
JAK2 (V617F) (JH1, JH2 Domain), GST-Tag, Avi-Tag, Biotin-Labeled (Dephosphorylated)	79115
JAK2 (V617F) (JH2 Domain), His-Avi-Tag, Biotin-Labeled	79075
JAK2 (V617F), FLAG-Avi-Tag, Biotin-Labeled (Dephosphorylated)	79426
JAK2 (V617F, W659A, W777A, F794H), His-Avi-Tag, Biotin-Labeled	79498
JAK2 (V617F, Y1007F, Y1008F)- (GSSG)8-EPOR	100709
JAK2 (W659A, W777A, F794H) (JH2 Domain), His-Avi-Tag, Biotin-Labeled	79463
JAK2, FLAG-Avi-Tag, Biotin-Labeled (Dephosphorylated)	79493
JAK3 (JH2 Domain), His-Tag	101270
JAK3, His-tag	40452
JNK1-β1(K55M), Kinase Deficient, GST-tag	40871
JNK2, GST-Tag	40113
KHS1 (GCKR), GST-tag	40115
LATS1, GST-Tag	79499
LATS2, GST-Tag	79042
LATS2, GST-Tag	100353
LCK, GST-tag	40470
LKB1/MO25/STRAD Complex	40069
Lok1 (Stk10), His-tag	40074
LRRK2 (D1994A), GST-Tag	40831
LRRK2 (G2019S), GST-Tag	40832
LRRK2 (G2385R), GST-Tag	40833
LRRK2 (I2020T), GST-Tag	40834
LRRK2 (R1441C), GST-Tag	40835
LRRK2 (R1441G), GST-Tag	40836
LRRK2 (R1441H), GST-Tag	40837
LRRK2 (Y1699C), GST-Tag	40838
LRRK2 (Y1699G), GST-Tag	40839
LRRK2, GST-Tag	40842
LTK, GST-Tag	101357
LYN A, GST-tag	40471
LYNB, GST-tag	40472
MAP4K3 (GLK), GST-tag	100624
MAPK10 (JNK3), GST-tag	40092
MAPK14 (p38α), His-Tag, Active	40243
MAPK14, Inactive, GST-tag	40091
MAPKAPK2, His-tag	40116
MAPKAPK2, Inactive, GST-Tag	40088

Proteins	Catalog#
MAPKAPK3, GST-tag	40117
MAPKAPK5, GST-tag	40118
MARK1, GST-tag	40119
MARK3, GST-tag	40120
MBP, His-Tag	40535
MEK1 (K97R), Inactive, His-tag	40075
MEK1, GST-tag	40123
MEK1, Non-activated, GST-tag	40531
MEK2, GST-tag	40125
MEKK2, GST-tag	40122
MEKK3, GST-tag	40124
MELK, GST-tag	40708
MINK1, GST-tag	40126
MKK4, His-Tag	101874
MKK6, GST-tag	40089
MLCK, GST-tag	40127
MLK1, GST-tag	41102
MLK2, GST-tag	41103
MLK3, GST-tag	41104
MLK4, GST-tag	41105
MNK1 (T385D), GST-tag	40078
MNK2, GST-tag	40128
Mouse AURORA A, GST-tag	40177
Mouse Cot (MAP3K8), His-GST-Tags	41050
Mouse EGFR, FLAG-Tag	40195
Mouse EPHB1, GST-tag	40199
Mouse FER, GST-tag	40419
Mouse ITK, GST-tag	40446
Mouse JNK1, GST-tag	40071
Mouse LC20, GST-tag	50127
Mouse MEK1 (EE), GST-tag	40121
Mouse PAK3, GST-tag	40147
Mouse PKM2, His-Tag	50296
Mouse Sphingosine kinase 1a, His-tag	40613
Mouse Sphingosine kinase 2, His-tag	40614
MRCKα, GST-tag	40129
MRCKβ, GST-tag	40130
MSK1, GST-Tag	40131

Proteins	Catalog#	Proteins	Catalog#
MSSK1 (SRPK3), GST-tag	40132	PDGFR $\alpha$ , GST-tag	40261
MST3, GST-tag	40133	PDGFR $\alpha$ , Mouse, GST-tag	40260
MST4, GST-tag	40134	PDGFR $\alpha$ (D842I), GST-Tag	100202
mTOR/Raptor/MLST8 Complex, FLAG-Tag, His-Tag (Human)	40300	PDGFR $\alpha$ (D842V), GST-Tag	79633
MUSK, His-tag	40079	PDGFR $\alpha$ (D842Y), GST-Tag	100201
MYLK2, His-tag	40135	PDGFR $\beta$ , His-tag	40263
MYLK3, GST-tag	40136	PDHK1, GST-Tag	40084
MYO3 $\beta$ , GST-tag	40137	PDK1, GST-tag	40080
NDR, GST-tag	40138	PDK3, GST-Tag	100927
NEK11, GST-tag	40139	PFKFB3, His-tag	71089
NEK2	40009	PHKG1, GST-tag	40150
NEK3, GST-tag	40140	PHKG2, GST-tag	40151
NEK6, His-tag	40014	PI3 Kinase (p110 $\alpha$ /p55 $\gamma$ ), His-Tag	40645
NEK7, GST-tag	40141	PI3 kinase (p110 $\alpha$ /p85 $\alpha$ )	40620
NEK7, His-Tag	100476	PI3 kinase (p110 $\alpha$ /p85 $\alpha$ )	40621
NEK9, GST-tag	40142	PI3 kinase (p110 $\beta$ /p85 $\alpha$ ), FLAG-tag	40622
Neuropilin-1, Avi-His-Tag	100870	PI3 kinase (p110 $\gamma$ /PIK3R5) complex, His-tag, GST-tag	40626
Neuropilin-1, Avi-His-Tag, Biotin-labeled, HiP™	100911	PI3 kinase (p110 $\delta$ /p85 $\alpha$ )	40628
NIK, GST-tag	40090	PI3 kinase (p120 $\gamma$ ), His-tag	40625
NUAK1, GST-Tag	101475	PI3 Kinase p110 $\alpha$ (E542K)/p85 $\alpha$ , FLAG-Tag	101913
NUAK2 (K81R), Inactive, GST-Tag	101474	PI3 Kinase p110 $\alpha$ (H1047L)/p85 $\alpha$ , FLAG-Tag	101914
NUAK2, GST-Tag	101476	PI3 kinase p110 $\alpha$ (N345K)/p85 $\alpha$	40646
p38 $\alpha$ , GST-tag	40070	PI3 kinase p110 $\alpha$ (E545K)/p85 $\alpha$	40640
p38 $\beta$ , GST-tag	40143	PI3 kinase p110 $\alpha$ (H1047R)/p85 $\alpha$ , FLAG-Tag	40641
p38 $\gamma$ (p38 $\gamma$ ), GST-tag	40144	PI3 kinase p110 $\alpha$ (K111N)/p85 $\alpha$	40639
p38 $\delta$ (p38 $\delta$ ), GST-tag	40145	PI3 Kinase [p110 $\alpha$ (E545K)/p85 $\alpha$ ], His-Tag	40644
p53 (Y220C), GST-Tag	101610	PIKFYVE, GST-Tag	101323
p53, GST-Tag	40511	PIM1, GST-tag	41107
p70S6K (S6K1), GST-tag	40062	PIM2, GST-tag	40153
p70S6Kb, GST-tag	40146	PIM3, Active, GST-tag (Mouse)	41108
p85 $\alpha$ , His-tag	40630	PIP5K1A, GST-tag	40305
PAK1	40072	PKA $\alpha$ , GST-tag	40154
PAK2	40073	PKA $\alpha$ , His-Tag	79044
PAK4	40076	PKA $\beta$ , GST-tag	40155
PAK5	40077	PKA $\gamma$ , GST-tag	40156
PAK6	40081	PKC $\iota/\lambda$ , GST-tag	40015
PAK7, GST-tag	40148	PKC $\alpha$ (PKCalpha), GST-tag	40157
PASK, GST-tag	40149	PKC $\beta$ I (PKCbeta1), GST-tag	40158

Proteins	Catalog#	Proteins	Catalog#
PKCβ II (PKCbeta2), GST-tag	40159	RIPK3, GST-Tag	79710
PKCγ (PKCgamma), GST-tag	40160	RIPK4, GST-Tag	79712
PKCδ (PKCdelta), GST-tag	40161	RIPK5, GST-Tag	79713
PKCε (PKCepsilon), GST-tag	40162	ROCK1, GST-tag	40085
PKCζ (PKCzeta), GST-tag	40163	ROCK2, GST-tag	40086
PKCη (PKCeta), GST-tag	40164	RON, GST-tag	40292
PKCθ (PKCtheta), GST-tag	40165	ROR1, Fc-Fusion (IgG1), Avi-Tag	79481
PKCμ (PKCmu), GST-tag	40166	ROR1, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79482
PKCν (PKCnu), GST-tag	40167	ROR1, Fc-Fusion (IgG1)-Avi-Tag, PE-labeled	100995
PKD2, GST-tag	40168	ROR1, GST-tag	40396
PKG1α, His-tag	40082	ROR2, Fc-Fusion (IgG1), Avi-Tag HiP™	100029
PKG1β, His-tag	40083	ROR2, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	100046
PKLR Var1 (PKR), His-tag	40500	ROR2, GST-tag	40296
PKLR Var2 (PKL), His-tag	40501	ROS1 (G2032R), GST-Tag	101640
PKM1 (PKM2/Variant 2), His-Tag	40502	ROS1 (G2101A), GST-Tag	101641
PKM2 (PKM2/Variant 1), His-Tag	50295	ROS1, GST-His-tag	40268
PKMYT1, GST-Tag	101569	RSK1, GST-tag	40172
PKN1 (PRK1), GST-Tag	79678	RSK2, GST-tag	40181
PKN2 (PRK2), GST-tag	40169	RSK3, GST-tag	40182
PKN3 (PRK3), GST-tag	101642	RSK4, GST-tag	40183
PLK1 Polo Box Domain, GST-Tag	50301	SGK1, GST-tag	40271
PLK1, His-tag	40033	SGK2, GST-tag	40272
PLK2, GST-tag	40034	SGK3, GST-tag	40273
PLK3 Polo Box Domain, MBP-tag, His-tag	50302	SIK (SNF1LK), His-tag	40274
PLK3, His-Tag	40035	SIK2, GST-tag	40264
PLK4, GST-tag	40236	SLK, GST-tag	40275
PRKG1, GST-tag	40175	SOK1, GST-tag	40276
PRKG2, GST-tag	40170	Sphingosine kinase 1, His-tag	40610
PRKX, GST-tag	40171	Sphingosine Kinase 2 (Transcript Variant 1), His-Tag	40612
PYK2, His-tag	40480	Sphingosine kinase 2, His-tag	40611
Rat ITK, GST-tag	40447	SRC, GST-tag	40483
Rat MET, GST-tag	40228	SRC, His-Tag	40484
RB, GST-tag	40595	SRPK1, GST-tag	40277
RBD-c-RAF, GST-Tag	100519	SRPK2, GST-tag	40278
Ret, GST-tag	40267	STK19, GST-Tag	100910
RIPK1, GST-Tag (Mouse)	100340	STK3(MST2), GST-tag	40013
RIPK1, GST-Th-Tag	40371	STK33, GST-tag	40093
RIPK2, GST-tag	40173	STK4(MST1), GST-tag	40017

Proteins	Catalog#	Proteins	Catalog#
STYXL1	30027	ULK2, GST-tag	40294
SYK, Catalytic, GST-Tag	40487	ULK3, His-tag	40295
SYK, Full length, GST-tag	40485	VEGFR2 (KDR), GST-Tag	40301
TAK1-TAB1, GST-tag	40279	VEGFR3 (FLT4), GST-tag	40302
TAOK1, GST-tag	40283	VPS34, FLAG-tag	40690
TAOK2, GST-tag	40284	VRK1, GST-tag	40397
TBK1, GST-Tag	40286	Wee1, FLAG-Tag	100154
TGFβR1, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled	100645	Wee1, GST-Th-Tag	40412
TGFβR1, Fc fusion, Avi-Tag	100644	WNK1, GST-Tag	100971
TGFβR1, GST-Tag	101575	WNK2, GST-Tag	100972
TGFβR2, Fc Fusion, Avi-tag (Mouse)	100676	WNK3, GST-Tag	100973
TGFβR2, Fc Fusion, Avi-tag	100628	WNK4, GST-Tag (Mouse)	100974
TGFβR2, Fc Fusion, Avi-tag, Biotin-labeled (Mouse)	100677	Yes1, GST-tag	40488
TGFβR2, GST-tag	40707	ZAK, GST-tag (Variant 2)	40297
Tie2, GST-tag	40370	ZAP70, GST-tag	40490
Tie2, His-tag	40270		
TLK2, GST-tag	40287	Substrates	Catalog#
TNIK, GST-Tag	11708	5X Raf Substrate	79569
TNK1, GST-Th-Tag	80019	CERK Substrate	40619
TOPK, GST-tag	40288	CERK Substrate	40618
TPM3-TRKA Mutant, GST-Tag	79080	IRS1-tide	79519
TRKA, GST-tag	40280	PI-3 Kinase Lipid Substrate	40560
TRKB, GST-tag	40281	Protein Tyrosine Kinase Substrate (poly-Glu,Tyr 4:1)	40217
TRKC (G623E) Mutant, Active)	40204	STAT5 Peptide	79864
TRKC (G623R L686M),Mutant, Active	40215		
TRKC (G623R) Mutant, Active)	40203		
TRKC (L686M), Mutant, Active	40216		
TRKC, GST-tag	40282		
TSSK1B, GST-tag	40289		
TSSK2, GST-tag	40290		
TTBK1, GST-Tag, His-Tag	40309		
TTK, GST-tag	40291		
TXK, GST-tag	40486		
TYK2 (JH2 domain), FLAG-Tag	100523		
TYK2, GST-Tag	100400		
Tyk2, His-Tag	40285		
TYRO3, GST-tag	40293		
ULK1, FLAG-Tag	40099		



## FIND A TRUSTED PARTNER FOR YOUR NEXT PROJECT



Ordering Information  
[orders@bpsbioscience.com](mailto:orders@bpsbioscience.com)



Technical Support  
[support@bpsbioscience.com](mailto:support@bpsbioscience.com)



US Sales Support  
[sales-team@bpsbioscience.com](mailto:sales-team@bpsbioscience.com)



International Sales Support  
[international@bpsbioscience.com](mailto:international@bpsbioscience.com)

BPS Bioscience, Inc.  
6405 Mira Mesa Blvd, Suite 100  
San Diego, CA 92121  
Tel: 858-202-1401

[bpsbioscience.com](http://bpsbioscience.com)

