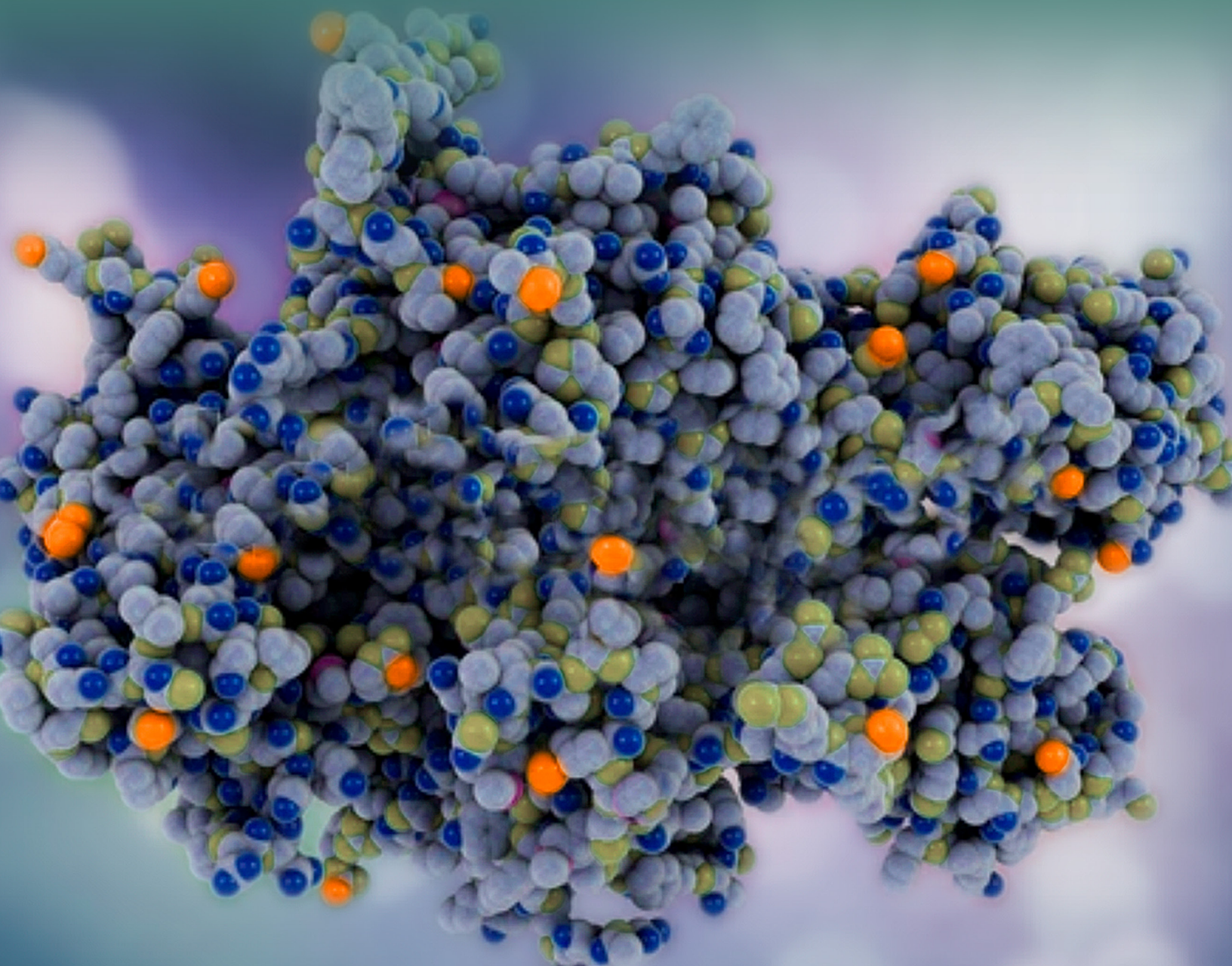




# PRODUCT SOLUTIONS FOR KINASES

Enabling Drug Discovery and Development Programs

Comprehensive Offering of Cellular Assays and Recombinant Enzymes  
for Drug Discovery Screening and Lead Optimization



# Comprehensive Kinase Solutions to Accelerate Your Kinase Drug Discovery

The Largest Commercial Offering of Optimized Kinase Enzymes and Cell-Based Assays for Drug Discovery Target Characterization, Screening and Profiling, and Lead Optimization Programs

Whether you are performing a biochemical enzymatic or functional cell-based assay screen, Eurofins DiscoverX is your drug discovery and development partner. As the new home of the original Millipore (Upstate) kinase enzyme portfolio covering over 70% of the human kinome combined with the clinically-proven PathHunter<sup>®</sup> mechanism-of-action (MOA)-reflective, cell-based assay platform, Eurofins DiscoverX has the kinase solutions you need to accelerate your drug discovery programs.



## RECOMBINANT KINASE ENZYMES HIGHLIGHTS

Access the most extensive commercial offering of high-quality recombinant active, inactive, and mutant kinases for screening and profiling potential drugs. Available in multiple pack sizes, including bulk sizing, to provide you with scalable options. Whether you want to perform all of your screening in-house or with KinaseProfiler services, you will receive the quality and confidence that comes with 100s of millions of data points generated over the past 30 years.

- **Comprehensive Portfolios** – Over 490 high quality active, inactive, and mutant kinase enzymes, including many exclusive enzymes such as the PIKK group (ATM, ATR, mTOR, and DNA-PK)
- **Optimal Expression Systems** – All recombinant kinases are produced in their optimal system (insect, mammalian, or bacterial) to give the correct characteristics for activity and inhibition
- **High Quality** – Rigorous quality control procedure ensures the highest purity (>95%), specific activity, and lot-to-lot consistency

## CELL-BASED KINASE ASSAY TYPES AND HIGHLIGHTS

MOA-reflective cell-based kinase assays for target characterization to lead optimization, and through clinical development to QC lot release.

### MOA-REFLECTIVE PathHunter CELL-BASED KINASE ASSAYS

Cell-based assays for analyzing ligand activation, receptor dimerization and internalization, SH2-recruitment, and screening of novel kinase inhibitors for receptor tyrosine kinases (RTK) and cytokine receptors. These assays provide cellular context to kinase activation to identify and profile novel therapeutic compounds and biologics.



- **Broadly Applicable** – Identify various ligands including anti-receptor, anti-ligand, or activating antibodies; non-ATP pocket binders (allosteric modulators); ligand binding inhibitors (ATP-competitors); or dimerization inhibitors
- **Accurate Reproducibility** – Superior quality, reproducible data with large assay windows and robust performance
- **East-to-Use** – Simple, homogeneous protocols with a chemiluminescent output that can be read on any benchtop luminometer.

## CELL-BASED KINASE ASSAY TYPES AND HIGHLIGHTS

### InCELL CELL-BASED TARGET ENGAGEMENT ASSAYS

InCELL Hunter™ and InCELL Pulse™ cellular compound-target engagement assays provide the ability to confirm compound cell entry and drug-target binding to intracellular targets or monitor drug-mediated protein degradation to assess compound efficacy and confirmation of MOA. The InCELL assay platform is ideal for screening inhibitors, validating hits identified in biochemical assays, measuring cellular EC<sub>50</sub> values, and ranking compounds in a native cellular environment.



- **Flexible** – Cellular, target engagement assays available as cell lines, ready-to-assay kits, or custom
- **Diverse** – Easily measure compound entry and drug-target binding, or accurately screen compounds in a high throughput format
- **Simple** – Binding assays that do not require custom chemical tracers, antibody reagents, or mass spec

### SPRINTER™ CELL-BASED TARGETED PROTEIN DEGRADATION ASSAYS

SPRINTER targeted protein degradation assays are for rapid screening of small molecule therapeutics and quantifying changes in endogenous protein levels in disease-relevant cell models. Detect protein turnover induced by targeted degrader molecules, such as PROTAC®s (Proteolysis Targeting Chimeras) or molecular glues, with higher sensitivity and more rapid kinetics than phenotypic endpoint assays (e.g. cell proliferation). Discover new molecular entities that modulate the endogenous levels of targeted proteins by using both SPRINTER platform with InCELL Pulse™ target engagement assays.



- **Homogeneous** – Simple, scalable, and homogeneous no-wash protocol amenable to high-throughput screening (HTS) formats for increased efficiency
- **Robust & Highly Sensitive** – Accurately detect target protein turnover at micro/nanomolar sensitivities
- **Rapid Results** – Obtain results in as little as 5 hours to select the right candidate and accelerate development programs
- **Optimized Signal** – EFC detection method features lysine-free tags to minimize artifacts

### ADP ACCUMULATION ASSAYS FOR KINASE PROFILING AND HTS APPLICATIONS

Homogeneous activity-based ADP accumulation assays are antibody-free screening; ideal for kinase profiling, high throughput screening (HTS), and identifying and characterizing phosphotransferase activity. In contrast to standard assays that rely on antibody detection of a phosphoepitope or monitoring ATP depletion as the result of kinase activity, these assays are gain-of-signal assays that generate a positive readout in direct proportion to ADP accumulation as a result of substrate phosphorylation activity.

- **Ideal Assays** – Kinetic and end-point modes ideal for inhibitor screening and profiling
- **Flexible** – Robust fluorescence readout allows for flexible assay read time
- **Simple** – Easy-to-use and universal for any phosphotransferases, such as ATPases, UTPases, and GTPases



## CUSTOM KINASE DEVELOPMENT CAPABILITIES

When an off-the-shelf option is unavailable, consider Eurofins DiscoverX custom cell lines, recombinant proteins, and assays optimized to your requirements.



- **Development Expertise** – Decades of cell-based assay development, cell line engineering, and recombinant enzyme development expertise
- **Cell Line Engineering Capability** – Exogenous expression approaches (constitutive vs inducible) or gene editing (e.g. KO/KI with CRISPR/Cas9)
- **Collaborative** – Consultative assay development with status updates through a dedicated project manager
- **Complete Solution** – Customized assay development with screening and profiling services within the same team

Learn more about Eurofins DiscoverX customer capabilities at [discoverx.com/CAD](https://discoverx.com/CAD)

---

## KINASE SCREENING AND PROFILING SERVICES

Screen your compounds against the broadest panel of wild-type and mutant human kinases to determine your compound potency and selectivity, and to advance your hits with the highest probability of being clinically effective and safe. Eurofins Discovery Services has over two decades of kinase screening and profiling experience with over 32,000 reports delivered screening for >1M kinase candidate compounds. With >1000 biochemical, including binding, activity, and cell-based kinase assays available, Eurofins Discovery Services partners with you to confidently confirm your hits.

- **Comprehensive** – Over 490 active-site directed competitive ligand binding assays (KINOMEscan® Technology) and 422 activity functional-based assays (KinaseProfiler® Technology) available to confirm activity assay hits
- **Conformational** – Confidently confirm your results across multiple biochemical and cellular orthogonal platforms
- **Flexible** – Customize panels specifically to your needs and modify as your program progresses

Learn more about kinase screening and profiling services at [eurofinsdiscoveryservices.com](https://eurofinsdiscoveryservices.com).

---

## PRODUCT TARGET LIST

The following tables indicates the available kinase **products**. For kinase services, please refer to our services website. The tables are organized by product type and includes all available kinase related toolbox products, biochemical kits, cell-based assays, and recombinant proteins. The cell-based assays section is organized by target and includes the product format (cell line, eXpress kit, bioassay kit, expression vector, and recombinant protein) indication. The recombinant proteins section is organized by target and includes product names (with reference of active, activated, inactive, or mutant state), expression system, tag, species, molecular weight, and catalog number.

### Eurofins Discovery Websites

**Products** [discoverx.com/kinases](https://discoverx.com/kinases)

**Services** [eurofinsdiscoveryservices.com](https://eurofinsdiscoveryservices.com)

Please refer to our websites for detailed information and the most update-to-date list of products currently available.

**Toolbox Products Parental Cell Lines, Starter Kit, and Expression Vectors**

Application	Expressed Protein	Catalog #	Product Name	Size	Target
Target Engagement Assay Development	Intracellular Targets	94-40075	InCELL Pulse™ Target Engagement Starter Kit	400 dp (4 X 96-well)	Kinases
Monitor SH2 Recruitment	mPLCG1(SH2)-EA	93-1124C3	PathHunter® UZOS mPLCG1(SH2)-EA Parental Cell Line	2 vials	RTKs
Monitor SH2 Recruitment	SHC1(SH2)-EA	93-1123C3	PathHunter® UZOS SHC1-EA Parental Cell Line	2 vials	RTKs
Translocation to the Membrane	MEM-EA	93-1101C3	PathHunter® UZOS MEM-EA Parental Cell Line	2 vials	RTKs
Translocation to the Endosome	ENDO-EA	93-1102C3	PathHunter® UZOS ENDO-EA Parental Cell Line	2 vials	RTKs
Cell-based Assay Development	Intracellular Targets	Varies	ProLabel™ & ProLink™ Expression Vectors	10 µg	RTKs, CTKs
Cell-based Assay Development	Intracellular Targets	Varies	InCELL Pulse™ Expression Vectors	10 µg	Kinases

**ADP Accumulation Assays**

Product Type	Target	Base Catalog #	Product Name	Size	Read-Out
Biochemical Kit	Any Kinase, ATPase,	90-0083	ADP Hunter™ Plus	Multiple	Fluorescent
Biochemical Kit	GTPase or UTPase	90-0071	ADP Quest™	Multiple	Fluorescent

**PathHunter® and InCELL Cell-Based Assays for RTKs, CTKs and Kinases**

*\* See last table for full recombinant protein list*

Application	Target	Kinase Type	Cell Line	eXpress Kit	Bioassay Kit	Expression Vector	Recombinant Protein*
Target Engagement	AAK1	Kinase				●	
Target Engagement	ABL1	Kinase		●		●	●
Target Engagement	ABL1(T315I)	Kinase				●	●
Target Engagement	ACVR1	Kinase				●	
Dimerization	ACVR1/ACVR2	RTK	●			●	
Dimerization	ACVR1B/BMPR2	RTK	●	●		●	
Dimerization	ACVR1C/ACVR2	RTK	●	●		●	
Dimerization	ACVR1C/ACVR2B	RTK	●	●		●	
Dimerization	ACVRL1/ACVR2	RTK	●	●		●	
Dimerization	ACVRL1/ACVR2B	RTK	●	●		●	
Dimerization	ACVRL1/BMPR2	RTK	●	●		●	
Target Engagement	AKT1	Kinase				●	●
Target Engagement	AURKA	Kinase				●	
Functional	AXL	RTK	●			●	●
Activity, Target Engagement	BLK	CTK	●	●		●	●
Dimerization	BMPR1A/ACVR2	RTK	●	●		●	
Dimerization	BMPR1A/ACVR2B	RTK	●			●	
Dimerization	BMPR1A/BMPR2	RTK	●	●		●	
Dimerization	BMPR1B/ACVR2A	RTK	●	●		●	
Dimerization	BMPR1B/ACVR2B	RTK	●			●	
Dimerization	BMPR1B/BMPR2	RTK	●			●	
Target Protein Degradation	BRAF(V600E)	Kinase	●	● <i>New</i>			
Target Engagement	BRAF	Kinase				●	●
Target Engagement	BTK	Kinase	●	●		●	●
Target Protein Degradation	BTK	Kinase	●	● <i>New</i>			
Target Engagement	BUB1	Kinase				●	
Target Engagement	CAMK2A	Kinase				●	●
Target Protein Degradation	CDKN1A (p21)	Kinase	●	● <i>New</i>			
Functional	cINSRa	RTK	●	●		●	
Functional	cINSRb	RTK	●	●		●	
Functional	c-KIT	RTK	●	●		●	●
Target Engagement	CLK1	Kinase	●	●		●	●
Target Engagement	CLK2	Kinase	●	●		●	●
Functional	c-MET	RTK				●	●
Dimerization	c-MET/c-MET	RTK				●	
Dimerization	c-MET/EGFR	RTK	● <i>New</i>			●	
Functional	c-Ret-GFRα1	RTK	●	●		●	●
Functional	c-Ret-GFRα2	RTK	●	●		●	●
Functional	c-Ret-GFRα3	RTK	●	●		●	
Target Engagement	CSF1R	Kinase	●	●		●	
Dimerization	CSF1R/CSF1R	RTK	●	●		●	
Dimerization	CSF2RB/CSF2RA	RTK	●	●		●	
Functional	CSF3R-JAK1	CTK	●	●		●	
Target Engagement	CSNK1D	Kinase	●	●		●	
Target Engagement	CSNK2A2	Kinase				●	
Functional	DDR1	RTK	●	●		●	●
Dimerization	EGFR/EGFR	RTK	●	●		●	
Dimerization	EGFR/ErbB2	RTK	●	●	●	●	
Dimerization	EGFR/ErbB3	RTK	●	●		●	
Functional	EphA4	RTK	●	●		●	●
Functional	EphA5	RTK	●	●		●	●
Functional	EphA7	RTK	●	●		●	●
Functional	EphB1	RTK	●	●		●	●
Functional	EphB2	RTK	●	●		●	●
Functional	EphB3	RTK	●	●		●	●
Functional	EphB4	RTK	●	●		●	●
Dimerization	EpoR/EpoR	RTK	●	●	●	●	
Functional	EpoR-JAK2	CTK	●	●		●	
Functional	ErbB1	RTK	●			●	

# Kinase Products Target List

PathHunter® and InCELL Cell-Based Assays for RTKs, CTKs and Kinases							
Application	Target	Kinase Type	Cell Line	eXpress Kit	Bioassay Kit	Expression Vector	Recombinant
Dimerization	ErbB2/ErbB3	RTK	●		●	●	
Functional	ErbB2-ErbB3	RTK	●	●		●	
Functional	ErbB4	RTK	●			●	●
Dimerization	ErbB4/ErbB4	RTK	●	●		●	
Target Engagement	ERK1	Kinase				●	
Target Engagement	FAK	Kinase				●	●
Functional	FGFR1	RTK	●	●		●	●
Functional	FGFR1- $\alpha$ -Klotho	RTK	●			●	
Functional	FGFR1- $\beta$ -Klotho	RTK	●			●	
Functional	FGFR1v3B	RTK	●			●	
Functional	FGFR2	RTK	●			●	●
Dimerization	FGFR3(G380R)/FGFR3(G380R)	RTK	●			●	
Dimerization	FGFR3/FGFR3	RTK	●			●	
Functional	FGFR4	RTK	●	●		●	●
Functional	FGFR4- $\alpha$ -Klotho	RTK	●			●	
Functional	FGFR4- $\beta$ -Klotho	RTK	●	●	●	●	
Activity, Target Engagement	FGR	CTK	●	●		●	●
Activity	Flt3	RTK	●	●		●	●
Activity	Flt4	RTK	●	●		●	●
Target Engagement	GAK	Kinase				●	
Functional	GHR-JAK1	CTK	●	●	●	●	
Functional	GHR-JAK2	CTK	●	●		●	
Target Engagement	HASPIN	Kinase				●	●
Target Engagement	HCK	Kinase	●	●		●	●
Target Engagement	HPK1	Kinase	●			●	
Functional, Target Engagement	IGF1R	RTK	●	●	●	●	
Functional	INSRa	RTK	●			●	
Functional	INSRb	RTK	●		●	●	
Functional	INSRa (mouse)	RTK	●	●		●	
Functional	INSRb (mouse)	RTK	●	●		●	
Target Engagement	JAK2(JH1)	Kinase				●	
Activity	JAK3	CTK	●	●		●	●
Activity	KDR	RTK	●	●		●	●
Dimerization	KDR/KDR	RTK	●	●	●	●	
Target Engagement	KIT	Kinase				●	●
Activity	LCK	CTK	●			●	●
Target Engagement	MEK1	Kinase	●	●		●	●
Target Engagement	p38 $\alpha$	Kinase				●	●
Target Engagement	PAK4	Kinase				●	●
Functional	PDGFR $\alpha$	RTK	●	●	●	●	●
Functional	PDGFR $\beta$	RTK	●	●	●	●	●
Target Engagement	PI3K $\delta$	Kinase	●			●	
Target Engagement	PIKFYVE	Kinase	●			●	
Target Engagement	PIM1	Kinase	●	●		●	
Target Engagement	PKAC $\alpha$	Kinase				●	
Target Engagement	PKC $\theta$	Kinase	●			●	
Target Engagement	PLK1	Kinase				●	●
Functional	PRLR-JAK1	CTK	●	●		●	
Functional	PRLR-JAK2	CTK	●	●		●	
Target Engagement	PYK2	Kinase	●	●		●	●
Target Engagement	RIPK1	Kinase	●			●	
Target Engagement	RIPK2	Kinase	●	●		●	●
Target Engagement	RIPK3	Kinase	●			●	
Target Engagement	SIK1	Kinase	●	●		●	●
Target Engagement	SRPK1	Kinase				●	●
Activity	SYK	CTK	●	●		●	●
Dimerization	TGFBR1/ACVR2	RTK	●	●		●	
Dimerization	TGFBR1/ACVR2B	RTK	●	●		●	
Dimerization	TGFBR1/TGFBR2	RTK	●	●		●	
Dimerization	TGFBR1/TGFBR2/ENG	RTK	●			●	
Functional	Tie2	RTK	●			●	●
Functional	TrkA	RTK	●	●		●	●
Functional	TrkA-P75	RTK	●	●	●	●	
Functional	TrkA-P75 (rat)	RTK	●	●		●	
Functional	TrkA-P75 (monkey)	RTK	●	●		●	
Functional	TrkB	RTK	●	●		●	●
Functional	TrkB-P75	RTK	●	●	●	●	
Functional	TrkC	RTK	●	●	●	●	●
Functional	TrkC-P75	RTK	●	●		●	
Activity	TYK2	CTK		●		●	●
Functional	TYRO3	RTK	●			●	
Target Engagement	ULK1	Kinase	● <i>New</i>			●	
Target Engagement	VPS34	Kinase				●	
Activity	YES1	CTK	●	●		●	●

Recombinant Kinases and Phosphatases		Multiple sizes including bulk sizes are available					
Product Name	Target	Target Type	Expression System	Tag	Species	Molecular Weight	Base Catalog No.
AAK1, active	AAK1	Kinase	Sf21 insect cells	His6	Human	43 kDa	16-043
Abl (H396P), active	Abl	Kinase	Sf21 insect cells	His6	Human	121 kDa	14-750
Abl (M351T), active	Abl	Kinase	Sf21 insect cells	His6	Human	121 kDa	14-757
Abl (Q252H), active	Abl	Kinase	Sf21 insect cells	His6	Human	121 kDa	14-751
Abl (T315I), active	Abl	Kinase	Sf21 insect cells	His6	Human	121 kDa	14-522
Abl (Y253F), active	Abl	Kinase	Sf21 insect cells	His6	Human	121 kDa	14-759
Abl, active	Abl	Kinase	Sf21 insect cells	His6	Human	121 kDa	14-529
Abl, active, mouse	Abl	Kinase	Sf21 insect cells	His6	Mouse	123.5 kDa	14-459
ACK1, active	ACK1	Kinase	Sf21 insect cells	GST	Human	71.5 kDa	14-756
ACTR2, active	ACTR2	Kinase	Sf21 insect cells	GST	Human	66 kDa	16-009
ActRIIB, active	ActRIIB	Kinase	Sf21 insect cells	GST	Human	61 kDa	16-052
Akt1/PKBα (6PH, S473D), active	Akt1/PKBα	Kinase	Sf21 insect cells	His6	Human	45 kDa	14-453
Akt1/PKBα, active	Akt1/PKBα	Kinase	Sf21 insect cells	His6	Human	59.9 kDa	14-276
Akt1/PKBα, unactive	Akt1/PKBα	Kinase	Sf21 insect cells	His6	Human	59 kDa	14-279
Akt2/PKBβ (6PH, S474D), active	Akt2/PKBβ	Kinase	Sf21 insect cells	His6	Human	42.8 kDa	14-339
Akt3/PKBγ (S472D), active	Akt3/PKBγ	Kinase	Sf21 insect cells	His6	Human	43 kDa	14-502
Alk, active	ALK	Kinase	Sf21 insect cells	His6	Human	63.8 kDa	14-555
ALK1, active	ALK1	Kinase	Sf21 insect cells	His6	Human	45 kDa	14-954
ALK2, active	ALK2	Kinase	Sf21 insect cells	His6	Human	45 kDa	14-937
ALK4, active	ALK4	Kinase	Sf21 insect cells	GST	Human	63.8 kDa	14-614
ALK6, active	ALK6	Kinase	Sf21 insect cells	His6	Human	44 kDa	14-941
AMPK (α1, β1, γ1), active	AMPK (α1, β1, γ1)	Kinase	E. coli	His6	Human	64 kDa (AMPKα1), 30 kDa (AMPKβ1), 38 kDa (AMPKγ1)	14-840
AMPK (α2, β1, γ1), active	AMPK (α2, β1, γ1)	Kinase	E. coli	His6	Human	63 kDa (AMPKα2), 30 kDa (AMPKβ1), 38 kDa (AMPKγ1)	14-902
A-Raf, active	A-Raf	Kinase	Sf21 insect cells	GST	Human	65 kDa	14-956
Arg, active	Arg	Kinase	Sf21 insect cells	His6	Human	121.9 kDa	14-521
Arg, active, mouse	Arg	Kinase	Sf21 insect cells	His6	Mouse	124.5 kDa	14-460
ARK5, active	NuaK1 (ARK5)	Kinase	Sf21 insect cells	His6	Human	78 kDa	14-661
ASK1, active	Ask1	Kinase	E. coli	GST, His6	Human	61.5 kDa	14-606
ATM, active	ATM	Kinase	Mammalian cell line	FLAG	Human	352 kDa	14-933
ATR/ATRIP, active	ATR/ATRIP	Kinase	Mammalian cell line	FLAG, cMyc	Human	303 kDa (ATR), 87 kDa (ATRIP)	14-953
Aurora A, active	Aurora-A	Kinase	Sf21 insect cells	His6	Human	46.9 kDa	14-511
Aurora B, active	Aurora-B	Kinase	Sf21 insect cells	GST	Human	40 kDa (Aurora B), 38 kDa (INCENP)	14-835
Aurora C, active	Aurora C	Kinase	Sf21 insect cells	GST, His6	Human	36 kDa (Aurora C), 38 kDa (INCENP)	14-911
Axl, active	Axl	Kinase	Sf21 insect cells	His6	Human	48 kDa	14-512
BIKe, active	BIKe	Kinase	Sf21 insect cells	His6	Human	44 kDa	16-044
Blk, active	Blk	Kinase	Sf21 insect cells	His6	Human	58.8 kDa	14-517
Blk, active, mouse	Blk	Kinase	Sf21 insect cells	His6	Mouse	57.7 kDa	14-316
BMPr2, active	BMPr2	Kinase	Sf21 insect cells	His6	Human	67 kDa	16-001
Bmx, active	Bmx	Kinase	Sf21 insect cells	His6	Human	79 kDa	14-499
B-Raf (V599E), active	B-Raf	Kinase	Sf21 insect cells	GST	Human	67.3 kDa	14-557
B-Raf (G1-415), active	B-Raf	Kinase	Sf21 insect cells	GST	Human	67.2 kDa	14-530
Brk, active	BRK	Kinase	Sf21 insect cells	His6	Human	55.3 kDa	14-613
BrSK1, active	BrSK1	Kinase	Sf21 insect cells	His6	Human	88.4 kDa	14-675
BrSK2, active	BrSK2	Kinase	Sf21 insect cells	His6	Human	78.5 kDa	14-655
BTK (E41K), active	BTK	Kinase	Sf21 insect cells	His6	Human	78.4 kDa	14-773
BTK (R28H), active	BTK	Kinase	Sf21 insect cells	His6	Human	78.4 kDa	14-765
BTK, active	BTK	Kinase	Sf21 insect cells	His6	Human	78.4 kDa	14-552
CaM Kinase I γ, active	CaM Kinase I γ	Kinase	Sf21 insect cells	GST	Human	80 kDa	14-967
CaM Kinase I, active	CaMKI	Kinase	E. coli	GST	Human	68 kDa	14-663
CaM Kinase Iβ, active	CaM Kinase Iβ	Kinase	Sf21 insect cells	GST	Human	65 kDa	15-001
CaM Kinase IIα, active	CaM Kinase IIα	Kinase	Sf21 insect cells	His6	Human	55 kDa	14-962
CaM Kinase IIβ, active	CaMKIIβ	Kinase	Sf21 insect cells	His6	Human	39.6 kDa	14-718
CaM Kinase IIγ, active	CaMKIIγ	Kinase	Sf21 insect cells	His6	Human	41.2 kDa	14-719
CaM Kinase IIδ, active	CaMKIIδ	Kinase	Sf21 insect cells	His6	Human	58 kDa	14-723
CaM Kinase IV, active	CaMKIV	Kinase	Sf21 insect cells	His6	Human	52.9 kDa	14-547
CaM Kinase Iδ, active	CaMKIδ	Kinase	Sf21 insect cells	His6	Human	46.7 kDa	14-731
CaMKK1, active	CaMKK1	Kinase	Sf21 insect cells	His6	Human	60 kDa	15-022
CaMKK2, active	CaMKK2	Kinase	Sf21 insect cells	His6	Human	64 kDa	14-931
Casein Kinase 1γ1, active	CK1γ1	Kinase	Sf21 insect cells	His6	Human	41.9 kDa	14-711
Casein Kinase 1γ2, active	CK1γ2	Kinase	Sf21 insect cells	His6	Human	49.7 kDa	14-712
Casein Kinase 1γ3, active	CK1γ3	Kinase	Sf21 insect cells	His6	Human	42.2 kDa	14-713
Casein Kinase 1δ (aa 1-294), active	CK1γ6	Kinase	E. coli	GST	Human	61 kDa	14-520
Casein Kinase 2, active	CK2	Kinase	Sf21 insect cells	GST, His6	Human	49 kDa (α-subunit), 53 kDa (β-subunit)	14-197
Casein Kinase 2α, active	CK2α	Kinase	Sf21 insect cells	His6	Human	48.7 kDa	14-445
Casein Kinase 2α2, active	CK2α2	Kinase	Sf21 insect cells	His6	Human	45.8 kDa	14-689
CD45	CD45	Phosphatase	E. coli	His6	Human	84 kDa	14-618
Cdc7/cyclin B1, active	Cdc7/cyclin B1	Kinase	Sf21 insect cells	GST	Human	92 kDa (Cdc7), 75 kDa (cyclin B1)	16-025
Cdk1/cyclin B, active	CDK1/cyclinB	Kinase	Sf21 insect cells	GST, His6	Human	35 kDa (cdk1), 75 kDa (cyclin B)	14-450
CDK12/cyclin K, active	CDK12/cyclin K	Kinase	Sf21 insect cells	FLAG/untagged	Human	165 kDa/65 kDa	16-008-K
CDK13/cyclin K, active	CDK13/cyclin K	Kinase	Sf21 insect cells	FLAG/untagged	Human	167 kDa/65 kDa	16-023-K
CDK14/cyclin Y, active	CDK14/cyclin Y	Kinase	Sf21 insect cells	His6/His6	Human	54 kDa/44 kDa	15-034-K
CDK16/cyclin Y, active	CDK16/cyclin Y	Kinase	Sf21 insect cells	GST/GST	Human	83 kDa/67 kDa	16-041
CDK17/cyclin Y, active	CDK17/cyclin Y	Kinase	Sf21 insect cells	GST/GST	Human	87 kDa/67 kDa	16-042
CDK18/cyclin Y, active	CDK18/cyclin Y	Kinase	Sf21 insect cells	GST/GST	Human	81 kDa/67 kDa	15-031
Cdk2/Cyclin A, active	CDK2/cyclinA	Kinase	Sf21 insect cells	GST, His6	Human	35 kDa	14-448
Cdk2/Cyclin E, active	CDK2/cyclinE	Kinase	Sf21 insect cells	GST, His6	Human	34 kDa (cdk2), 74 kDa (cyclin)	14-475
Cdk3/Cyclin E, active	CDK3/cyclinE	Kinase	Sf21 insect cells	GST, His6	Human	36 kDa (cdk3), 74 kDa (cyclin E)	14-487
CDK4/cyclin D3, active	CDK4/cyclin D3	Kinase	Sf21 insect cells	GST	Human	61 kDa (CDK4), 59 kDa (cyclin D3)	14-957
Cdk5/p25, active	CDK5/p25	Kinase	Sf21 insect cells	GST, His6	Human	34.4 kDa (cdk5), 49.4 kDa (p25)	14-516

Recombinant Kinases and Phosphatases								Multiple sizes including bulk sizes are available
Product Name	Target	Target Type	Expression System	Tag	Species	Molecular Weight	Base Catalog No.	
Cdk5/p35, active	CDK5/p35	Kinase	Sf21 insect cells	GST, His6	Human	34 kDa (cdk5), 61 kDa (p35 with GST-tag)	14-477	
Cdk6/Cyclin D3, active	CDK6/cyclinD3	Kinase	Sf21 insect cells	GST, His6	Human	38 kDa (cdk6), 59 kDa (cyclin D3)	14-519	
Cdk7/Cyclin H/MAT1 (CAK complex), active	CDK7/cyclinH/MAT1	Kinase	Sf21 insect cells	GST, His6	Human	39.9 kDa (cdk7), 37.8 kDa (cyclin H), 62.8 kDa (MAT1)	14-476	
Cdk9/Cyclin T1, active	CDK9/cyclin T1	Kinase	Sf21 insect cells	His6	Human	44 kDa (cdk9), 80.79 kDa (cyclinT1)	14-685	
CDKL1, active	CDKL1	Kinase	Sf21 insect cells	His6	Human	39 kDa	16-002-K	
CDKL2, active	CDKL2	Kinase	Sf21 insect cells	His6	Human	60 kDa	16-034	
CDKL3, active	CDKL3	Kinase	Sf21 insect cells	His6	Human	55 kDa	16-027	
CDKL4, active	CDKL4	Kinase	Sf21 insect cells	His6	Human	40 kDa	16-031	
ChaK1, active	ChaK1	Kinase	Sf21 insect cells	His6	Human	82 kDa	14-961	
CHK1, active	CHK1	Kinase	Sf21 insect cells	GST	Human	81.6 kDa	14-346	
CHK2 (hu,5-end, R145W)	CHK2	Kinase	E. coli	GST, His6	Human	89.6 kDa	14-740	
CHK2 (i157T), active	CHK2	Kinase	E. coli	GST, His6	Human	89.6 kDa	14-741	
CHK2, active	CHK2	Kinase	E. coli	GST, His6	Human	89.6 kDa	14-347	
CK1 $\alpha$ , active	CK1 $\alpha$	Kinase	Sf21 insect cells	FLAG, His6	Human	42 kDa	16-050-K	
CK1 $\epsilon$ , active	CK1 $\epsilon$	Kinase	Sf21 insect cells	GST	Human	67 kDa	16-045-K	
c-Kit (D816H), active	c-Kit	Kinase	Sf21 insect cells	GST	Human	76.7 kDa	14-726	
c-Kit (D816V), active	c-Kit	Kinase	Sf21 insect cells	GST	Human	76.7 kDa	14-611	
c-Kit (V560G), active	c-Kit	Kinase	Sf21 insect cells	GST	Human	76.6 kDa	14-730	
c-Kit (V654A), active	c-Kit	Kinase	Sf21 insect cells	GST	Human	76.7 kDa	14-733	
c-Kit, active	c-Kit	Kinase	Sf21 insect cells	GST	Human	76.7 kDa	14-559	
CLK1, active	CLK1	Kinase	Sf21 insect cells	His6	Human	45 kDa	15-026	
CLK1, active	CLK1	Kinase	Sf21 insect cells	His6	Human	45 kDa	14-920	
CLK2, active	CLK2	Kinase	Sf21 insect cells	GST	Human	70.4 kDa	14-774	
CLK3, active	CLK3	Kinase	Sf21 insect cells	His6	Human	63 kDa	14-724	
CLK4, active	CLK4	Kinase	Sf21 insect cells	His6	Human	46 kDa	14-917	
CRIK, active	CRIK	Kinase	Sf21 insect cells	GST	Human	77 kDa	16-040	
CSK, active	CSK	Kinase	E. coli	GST	Human	77.6 kDa	14-458	
DAPK1, active	DAPK1	Kinase	Sf21 insect cells	GST	Human	60.9 kDa	14-692	
DAPK2, active	DAPK2	Kinase	Sf21 insect cells	GST	Human	70.6 kDa	14-657	
DCAMKL1, active	DCAMKL1	Kinase	Sf21 insect cells	His6	Human	85 kDa	14-986	
DCAMKL2, active	DCAMKL2	Kinase	Sf21 insect cells	His6	Human	39.1 kDa	14-716	
DCAMKL3, active	DCAMKL3	Kinase	Sf21 insect cells	His6	Human	38 kDa	14-943	
DDR1, active	DDR1	Kinase	Sf21 insect cells	His6	Human	50 kDa	14-942	
DDR2, active	DDR2	Kinase	Sf21 insect cells	His6	Human	48 kDa	14-579	
DMPK, active	DMPK	Kinase	Sf21 insect cells	His6	Human	65 kDa	14-649	
DNA-PK, active	DNA-PK	Kinase	Mammalian cell line	FLAG	Human	470 kDa	14-950	
DRAK1, active	DRAK1	Kinase	Sf21 insect cells	His6	Human	50.4 kDa	14-668	
DRAK2, active	DRAK2	Kinase	Sf21 insect cells	His6	Human	46 kDa	16-003	
DUSP22	DUSP22	Phosphatase	E. coli	GST	Human	47.3 kDa	14-641	
DYRK1A, active	DYRK1A	Kinase	Sf21 insect cells	GST	Human	112 kDa	14-951	
DYRK1B, active	DYRK1B	Kinase	Sf21 insect cells	GST	Human	96 kDa	14-944	
DYRK2, active	DYRK2	Kinase	Sf21 insect cells	His6	Human	63.5 kDa	14-669	
DYRK3, active	DYRK3	Kinase	Sf21 insect cells	GST	Human	94 kDa	15-002	
eEF-2K, active	eEF-2K	Kinase	E. coli	GST	Human	108.5 kDa	14-654	
EGFR (L858R), active	EGFR	Kinase	Sf21 insect cells	GST	Human	85.8 kDa	14-626	
EGFR (L861Q), active	EGFR	Kinase	Sf21 insect cells	GST	Human	85.8 kDa	14-627	
EGFR (T790M), active	EGFR	Kinase	Sf21 insect cells	GST	Human	86 kDa	14-725	
EGFR (T790M, L858R), active	EGFR	Kinase	Sf21 insect cells	GST	Human	85.8 kDa	14-721	
EGFR, active	EGFR	Kinase	Sf21 insect cells	GST	Human	86 kDa	14-531	
EphA1, active	EphA1	Kinase	Sf21 insect cells	His6	Human	50.6 kDa	14-653	
EphA2, active	EphA2	Kinase	Sf21 insect cells	His6	Human	38 kDa	14-560	
EphA3, active	EphA3	Kinase	Sf21 insect cells	His6	Human	49 kDa	14-644	
EphA4, active	EphA4	Kinase	Sf21 insect cells	His6	Human	37 kDa	14-574	
EphA5, active	EphA5	Kinase	Sf21 insect cells	His6	Human	37.9 kDa	14-639	
EphA7, active	EphA7	Kinase	Sf21 insect cells	His6	Human	38 kDa	14-672	
EphA8, active	EphA8	Kinase	Sf21 insect cells	His6	Human	37.1 kDa	14-673	
EphB1, active	EphB1	Kinase	Sf21 insect cells	His6	Human	51.4 kDa	14-674	
EphB2, active	EphB2	Kinase	Sf21 insect cells	His6	Human	52 kDa	14-553	
EphB3, active	EphB3	Kinase	Sf21 insect cells	His6	Human	40 kDa	14-561	
EphB4, active	EphB4	Kinase	Sf21 insect cells	His6	Human	51 kDa	14-554	
ErbB2, active	ErbB2	Kinase	Sf21 insect cells	His6	Human	69 kDa	14-939	
ErbB4, active	ErbB4	Kinase	Sf21 insect cells	His6	Human	36.1 kDa	14-569	
FAK, active	FAK	Kinase	Sf21 insect cells	His6	Human	35.5 kDa	14-720	
Fer, active	Fer	Kinase	Sf21 insect cells	His6	Human	35.6 kDa	14-605	
Fes/Fps, active	Fes/Fps	Kinase	Sf21 insect cells	His6	Human	98 kDa	14-473	
FGFR1 (V561M), active	FGFR1	Kinase	Sf21 insect cells	GST	Human	62.3 kDa	14-734	
FGFR1, active	FGFR1	Kinase	Sf21 insect cells	GST	Human	62 kDa	14-582	
FGFR2 (N549H), active	FGFR2	Kinase	Sf21 insect cells	His6	Human	38.1 kDa	14-742	
FGFR2, active	FGFR2	Kinase	Sf21 insect cells	His6	Human	38 kDa	14-617	
FGFR3, active	FGFR3	Kinase	Sf21 insect cells	His6	Human	36.8 kDa	14-464	
FGFR4, active	FGFR4	Kinase	Sf21 insect cells	His6	Human	36.4 kDa	14-583	
Fgr, active	Fgr	Kinase	Sf21 insect cells	His6	Human	60.4 kDa	14-568	
Flt1, active	Flt1	Kinase	Sf21 insect cells	GST	Human	89 kDa	14-923	
Flt-3 (D835Y), active	Flt3	Kinase	Sf21 insect cells	GST	Human	77.4 kDa	14-610	
Flt-3, active	Flt3	Kinase	Sf21 insect cells	GST	Human	77.4 kDa	14-500	
Flt-4, active	Flt4	Kinase	Sf21 insect cells	GST	Human	90.9 kDa	14-681	
Fms (Y969C), active	Fms	Kinase	Sf21 insect cells	His6	Human	50.2 kDa	14-820	
Fms, active	Fms	Kinase	Sf21 insect cells	His6	Human	50.2 kDa	14-551	
Fyn, active	Fyn	Kinase	Sf21 insect cells	His6	Human	61.8 kDa	14-441	
GAK, active	GAK	Kinase	Sf21 insect cells	FLAG, His6	Human	40.2 kDa	16-055	
GCK, active	GCK	Kinase	Sf21 insect cells	GST	Human	80.6 kDa	14-743	
GCN2, active	GCN2	Kinase	Sf21 insect cells	GST	Human	215 kDa	14-934	
GRK1, active	GRK1	Kinase	Sf21 insect cells	His6	Human	68 kDa	14-935	
GRK2, active	RK2	Kinase	Sf21 insect cells	His6	Human	82 kDa	14-965	
GRK3, active	GRK3	Kinase	Sf21 insect cells	His6	Human	81 kDa	15-028	
GRK4, active	GRK4	Kinase	Sf21 insect cells	His6	Human	67.7 kDa	16-057	



# Kinase & Phosphatase Products Target List

Email [customerservicedrx@eurofins.com](mailto:customerservicedrx@eurofins.com) to request custom products ending in Catalog No. -K

2024

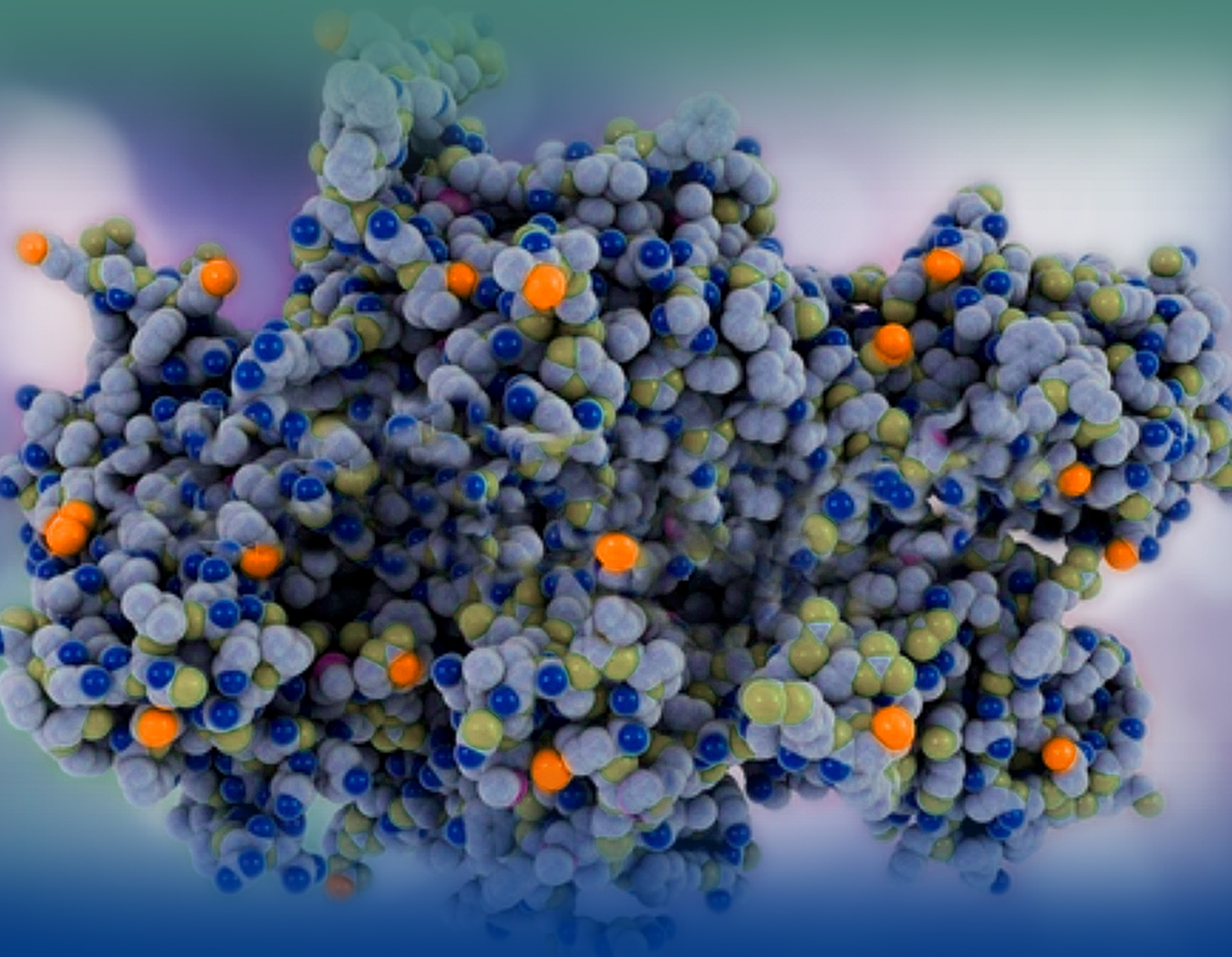
Recombinant Kinases and Phosphatases		Multiple sizes including bulk sizes are available					
Product Name	Target	Target Type	Expression System	Tag	Species	Molecular Weight	Base Catalog No.
GRK5, active	GRK5	Kinase	Sf21 insect cells	His6	Human	71.6 kDa	14-714
GRK6, active	GRK6	Kinase	Sf21 insect cells	His6	Human	69.8 kDa	14-715
GRK7, active	GRK7	Kinase	Sf21 insect cells	His6	Human	66 kDa	14-752
GSK3α, active	GSK3α	Kinase	Sf21 insect cells	His6	Human	55 kDa	14-492
GSK3β, active	GSK3β	Kinase	Sf21 insect cells	His6	Human	51 kDa	14-306
Haspin, active	haspin	Kinase	Sf21 insect cells	His6	Human	41.1 kDa	14-744
Hck, activated	Hck	Kinase	Sf21 insect cells	His6	Human	34.1 kDa	14-843
Hck, active	Hck	Kinase	Sf21 insect cells	His6	Human	34.1 kDa	14-577
HePTP	HePTP	Phosphatase	E. coli	His6	Human	40.7 kDa	14-593
HIPK1, active	HIPK1	Kinase	Sf21 insect cells	His6	Human	49.6 kDa	14-679
HIPK2, active	HIPK2	Kinase	Sf21 insect cells	His6	Human	49.6 kDa	14-623
HIPK3, active	HIPK3	Kinase	Sf21 insect cells	His6	Human	49.7 kDa	14-680
HIPK4, active	HIPK4	Kinase	Sf21 insect cells	His6	Human	71 kDa	15-003
HPK1, active	HPK1	Kinase	Sf21 insect cells	GST, His6	Human	67 kDa	14-968
HRI, active	HRI	Kinase	Sf21 insect cells	GST	Human	83 kDa	16-013
Histone H3	Histone	Substrate	E. coli	His6	Human	17.5 kDa	12-800
ICK, active	ICK	Kinase	Sf21 insect cells	His6	Human	40 kDa	15-023
IGF-1R, activated	IGF-1R	Kinase	Sf21 insect cells	His6	Human	48 kDa	14-802
IGF-1R (δ1-958), active	IGF-1R	Kinase	Sf21 insect cells	His6	Human	48 kDa	14-465
IKKα, active	IKKα	Kinase	Sf21 insect cells	GST	Human	111.6 kDa	14-461
IKKβ, active	IKKβ	Kinase	Sf21 insect cells	His6	Human	88 kDa	14-485
IKKe, active	IKKe	Kinase	Sf21 insect cells	GST, FLAG	Human	109 kDa	14-926
Insulin Receptor, activated	IR	Kinase	Sf21 insect cells	His6	Human	38.6 kDa	14-803
Insulin Receptor, active	IR	Kinase	Sf21 insect cells	His6	Human	36.8 kDa	14-466
IRAK1, active	IRAK1	Kinase	Sf21 insect cells	His6	Human	59.7 kDa	14-684
IRAK4, active	IRAK4	Kinase	Sf21 insect cells	His6	Human	57.1 kDa	14-599
IRE1, active	IRE1	Kinase	Sf21 insect cells	His6	Human	62 kDa	14-930
IRE2, active	IRE2	Kinase	Sf21 insect cells	His6	Human	56 kDa	16-054-K
IRR, active	IRR	Kinase	Sf21 insect cells	His6	Human	37.9 kDa	14-645
Itk, active	Itk	Kinase	Sf21 insect cells	His6	Human	34 kDa	14-660
JAK1, active	JAK1	Kinase	Sf21 insect cells	GST	Human	61 kDa	14-918
JAK2, active	JAK2	Kinase	Sf21 insect cells	His6	Human	38.9 kDa	14-640
JAK3, active	JAK3	Kinase	Sf21 insect cells	His6	Human	39.9 kDa	14-629
JNK1α1/SAPK1c, active	JNK1α1/SAPK1c	Kinase	Sf21 insect cells	His6	Human	45 kDa	14-327
JNK1α1/SAPK1c, unactive	JNK1α1/SAPK1c	Kinase	Sf21 insect cells	His6	Human	45 kDa	14-328
JNK2α2/SAPK1a, active	JNK2α2/SAPK1a	Kinase	Sf21 insect cells	His6	Human	49.2 kDa	14-329
JNK2α2/SAPK1a, unactive	JNK2α2/SAPK1a	Kinase	Sf21 insect cells	His6	Human	49.2 kDa	14-330
JNK3/SAPK1b, active	JNK3/SAPK1b	Kinase	Sf21 insect cells	His6	Human	53 kDa	14-501
JNK3/SAPK1b, unactive	JNK3/SAPK1b	Kinase	Sf21 insect cells	His6	Human	53 kDa	14-523
KDR, active	KDR	Kinase	Sf21 insect cells	His6	Human	67.9 kDa	14-630
Lambda PP, active, bacteriophage	Lambda PP	Phosphatase	E. coli	His6	Bacteriophage	26 kDa	14-946MG
LATS1, active	LATS1	Kinase	Sf21 insect cells	GST, His6	Human	90 kDa (LATS1), 30 kDa (MOBK1A)	14-988
LATS2, active	LATS2	Kinase	Sf21 insect cells	GST, His6	Human	89 kDa (LATS2), 30 kDa (MOBK1A)	14-987
Lck, activated	Lck	Kinase	Sf21 insect cells	His6	Human	59 kDa	14-842
Lck, active	Lck	Kinase	Sf21 insect cells	His6	Human	59 kDa	14-442
LIM Kinase 1, active	LIMK1	Kinase	Sf21 insect cells	His6	Human	41.2 kDa	14-656
LIM Kinase 1, unactive	LIMK1	Kinase	Sf21 insect cells	His6	Human	41.8 kDa	14-659
LIMK2, active	LIMK2	Kinase	Sf21 insect cells	His6	Human	76 kDa	16-032
LKB1/STRADA/MO25α, active	LKB1	Kinase	Sf21 insect cells	GST, His6	Human	51 kDa (LKB1), 76 kDa (STRADA), 66 kDa (MO25α)	14-596
LMPTP-A	LMPTP-A	Phosphatase	E. coli	GST	Human	44.5 kDa	14-619
LMPTP-B	LMPTP-B	Phosphatase	E. coli	GST	Human	44 kDa	14-620
LOK, active	LOK	Kinase	Sf21 insect cells	His6	Human	43.2 kDa	14-686
LRRK2, active	LRRK2	Kinase	Sf21 insect cells	FLAG	Human	180 kDa	14-919
LTK, active	LTK	Kinase	Sf21 insect cells	GST	Human	74 kDa	14-958
Lyn, active	Lyn	Kinase	Sf21 insect cells	His6	Human	59.6 kDa	14-510
Lyn, active, mouse	Lyn	Kinase	Sf21 insect cells	His6	Mouse	60 kDa	14-315
MAK, active	MAK	Kinase	Sf21 insect cells	GST	Human	80 kDa	16-010
MAP Kinase 1/Erk1, active	MAPK1	Kinase	E. coli	GST	Human	69.9 kDa	14-439
MAP Kinase 1/Erk1, unactive	MAPK1	Kinase	E. coli	GST	Human	70 kDa	14-515
MAP Kinase 2/Erk2, active	MAPK2	Kinase	E. coli	GST	Human	67.8 kDa	14-550
MAP Kinase 2/Erk2, active, mouse	MAPK2	Kinase	E. coli	GST	Mouse	67.8 kDa	14-173
MAP Kinase 2/Erk2, unactive	MAPK2	Kinase	E. coli	GST	Human	67.8 kDa	14-536
MAP Kinase 2/Erk2, unactive, mouse	MAPK2	Kinase	E. coli	GST	Mouse	67.8 kDa	14-198
MAP4K3, active	MAP4K3	Kinase	Sf21 insect cells	GST, His6	Human	64 kDa	15-030-K
MAP4K5, active	MAP4K5	Kinase	Sf21 insect cells	GST	Human	122 kDa	15-004-K
MAPK4, active	MAPK4	Kinase	Sf21 insect cells	His6	Human	66 kDa	15-012
MAPKAP Kinase 2, active	MAPKAP-K2	Kinase	E. coli	GST	Human	70.2 kDa	14-337
MAPKAP Kinase 2, unactive	MAPKAP-K2	Kinase	E. coli	GST	Human	66 kDa	14-349
MAPKAP Kinase 3, active	MAPKAP-K3	Kinase	E. coli	GST	Human	69.8 kDa	14-585
MAPKAP Kinase 3, unactive	MAPKAP-K3	Kinase	E. coli	GST	Human	69.8 kDa	14-586
MARK1, active	MARK1	Kinase	Sf21 insect cells	His6	Human	92.7 kDa	14-651
MARK3, active	MARK3	Kinase	Sf21 insect cells	GST	Human	108 kDa	15-032-K
MARK4, active	MARK4	Kinase	Sf21 insect cells	GST	Human	109 kDa	15-009-K
MATK, active	MATK	Kinase	Sf21 insect cells	His6	Human	53 kDa	16-056
MEK1, active	MEK1	Kinase	E. coli	GST, His6	Human	71 kDa	14-429
MEK1, unactive	MEK1	Kinase	E. coli	GST, His6	Human	71 kDa	14-420
MEK2, active	MEK2	Kinase	Sf21 insect cells	His6	Human	46 kDa	14-528
MEK2, unactive	MEK2	Kinase	Sf21 insect cells	His6	Human	46 kDa	14-532
MEKK2, active	MEKK2	Kinase	Sf21 insect cells	His6	Human	74 kDa	14-963
MEKK3, active	MEKK3	Kinase	Sf21 insect cells	GST	Human	98 kDa	16-015-K
MELK, active	MELK	Kinase	Sf21 insect cells	GST	Human	66 kDa	14-909
Mer, active	Mer	Kinase	Sf21 insect cells	GST	Human	64.4 kDa	14-728
Met (D1246H), active	Met (D1246H)	Kinase	Sf21 insect cells	His6	Human	50.1 kDa	14-915
Met (D1246N), active	Met	Kinase	Sf21 insect cells	His6	Human	50.1 kDa	14-818
Met (M1268T), active	Met	Kinase	Sf21 insect cells	His6	Human	50.1 kDa	14-817
Met (Y1248C), active	Met	Kinase	Sf21 insect cells	His6	Human	50 kDa	14-805
Met (Y1248D), active	Met	Kinase	Sf21 insect cells	His6	Human	50.0 kDa	14-816
Met (Y1248H), active	Met	Kinase	Sf21 insect cells	His6	Human	50.1 kDa	14-804

Recombinant Kinases and Phosphatases		Multiple sizes including bulk sizes are available					
Product Name	Target	Target Type	Expression System	Tag	Species	Molecular Weight	Base Catalog No.
Met, active	Met	Kinase	Sf21 insect cells	His6	Human	50 kDa	14-526
MINK, active	MINK	Kinase	Sf21 insect cells	His6	Human	38.7 kDa	14-615
MKK3, active	MKK3	Kinase	Sf21 insect cells	His6	Human	37 kDa	16-037-K
MKK4/SKK1, active, mouse	MKK4	Kinase	E. coli	GST	Mouse	67.7 kDa	14-377
MKK4/SKK1, inactive, mouse	MKK4	Kinase	E. coli	GST	Mouse	65 kDa	14-378
MKK6/SKK3 (S599D, T603D), active	MKK6	Kinase	E. coli	MBP	Human	80.6 kDa	14-537
MKK6/SKK3, active	MKK6	Kinase	E. coli	MBP	Human	82 kDa	14-303
MKK6/SKK3, inactive	MKK6	Kinase	E. coli	Mal-E	Human	80.6 kDa	14-304
MKP5, active	MKP5	Phosphatase	E. coli	His6	Human	20.6 kDa	14-779
MLCK, active	MLCK	Kinase	Sf21 insect cells	His6	Human	43.7 kDa	14-638
MLK1, active	MLK1	Kinase	Sf21 insect cells	His6	Human	35.3 kDa	14-690
MLK2, active	MLK2	Kinase	Sf21 insect cells	His6	Human	54 kDa	14-964
MLK3, active	MLK3	Kinase	Sf21 insect cells	GST	Human	79 kDa	16-039-K
MLK4, active	MLK4	Kinase	Sf21 insect cells	GST	Human	61 kDa	16-051
Mnk2, active	Mnk2	Kinase	Sf21 insect cells	His6	Human	50 kDa	14-664
MOK, active	MOK	Kinase	Sf21 insect cells	GST, His6	Human	70 kDa	14-960
MRCK $\alpha$ , active	MRCK $\alpha$	Kinase	Sf21 insect cells	FLAG, His6	Human	56 kDa	16-046-K
MRCK $\alpha$ , active	MRCK $\alpha$	Kinase	Sf21 insect cells	His6	Human	55.3 kDa	14-691
MRCK $\beta$ , active	MRCK $\beta$	Kinase	Sf21 insect cells	His6	Human	55.5 kDa	14-643
MSK1, active	MSK1	Kinase	Sf21 insect cells	His6	Human	94 kDa	14-548
MSK1, active (flag tag)	MSK1	Kinase	Sf21 insect cells	His6, FLAG	Human	94 kDa	14-438
MSK2, active	MSK2	Kinase	Sf21 insect cells	His6	Human	89.9 kDa	14-616
MSK2, inactive	MSK2	Kinase	Sf21 insect cells	His6	Human	89.9 kDa	14-625
MSSK1, active	MSSK1	Kinase	Sf21 insect cells	His6	Human	65.7 kDa	14-665
MST1, active	MST1	Kinase	Sf21 insect cells	His6	Human	57 kDa	14-624
MST2, active	MST2	Kinase	Sf21 insect cells	His6	Human	57.2 kDa	14-524
MST3, active	MST3	Kinase	Sf21 insect cells	GST	Human	61.2 kDa	14-695
MST4, active	MST4	Kinase	Sf21 insect cells	His6	Human	38 kDa	14-928
mTOR (I362-end), active	mTOR (FRAP1)	Kinase	Sf21 insect cells	FLAG	Human	137.3 kDa	14-770
MuSK, active	MuSK	Kinase	Sf21 insect cells	His6	Human	43.1 kDa	14-634
MYLK2, active	MYLK2	Kinase	Sf21 insect cells	His6	Human	66 kDa	14-966
MYO3B, active	MYO3B	Kinase	Sf21 insect cells	GST	Human	65 kDa	15-008-K
NDR1, active	NDR1	Kinase	Sf21 insect cells	GST, His6	Human	83 kDa (NDR1), 30 kDa (MOBK11A)	14-989
NDR2, active	NDR2	Kinase	Sf21 insect cells	GST, His6	Human	72kDa	16-028-K
NEK1, active	NEK1	Kinase	Sf21 insect cells	GST, His6	Human	86 kDa	15-020
NEK11, active	NEK11	Kinase	Sf21 insect cells	His6	Human	59.3 kDa	14-700
NEK2, active	NEK2	Kinase	Sf21 insect cells	His6	Human	52.8 kDa	14-545
NEK3, active	NEK3	Kinase	Sf21 insect cells	His6	Human	62.2 kDa	14-694
NEK4, active	NEK4	Kinase	Sf21 insect cells	His6	Human	96 kDa	15-033-K
NEK5, active	NEK5	Kinase	Sf21 insect cells	His6	Human	39.9 kDa	15-037
NEK6, active	NEK6	Kinase	Sf21 insect cells	His6	Human	40.5 kDa	14-578
NEK7, active	NEK7	Kinase	Sf21 insect cells	His6	Human	37.9 kDa	14-565
NEK9, active	NEK9	Kinase	Sf21 insect cells	His6	Human	41 kDa	14-936
NIM1, active	NIM1	Kinase	Sf21 insect cells	GST	Human	77 kDa	14-945
NLK, active	NLK	Kinase	Sf21 insect cells	His6	Human	48.7 kDa	14-676
NUAK2, active	NUAK2	Kinase	Sf21 insect cells	GST	Human	96 kDa	15-005
OSR1, active	OSR1	Kinase	Sf21 insect cells	GST	Human	86 kDa (OSR1), 67 kDa (MO25 $\alpha$ )	16-048
p38 $\alpha$ /SAPK2 $\alpha$ (T106M), active	p38 $\alpha$ /SAPK2 $\alpha$	Kinase	E. coli	GST	Human	67.7 kDa	14-687
p38 $\alpha$ /SAPK2 $\alpha$ , active	p38 $\alpha$ /SAPK2 $\alpha$	Kinase	E. coli	GST	Human	67.7 kDa	14-251
p38 $\alpha$ /SAPK2 $\alpha$ , inactive	p38 $\alpha$ /SAPK2 $\alpha$	Kinase	E. coli	GST	Human	67.7 kDa	14-252
p38 $\beta$ 2/SAPK2b2, active	p38 $\beta$ 2/SAPK2b2	Kinase	E. coli	GST	Human	71 kDa	14-253
p38 $\beta$ 2/SAPK2b2, inactive	p38 $\beta$ 2/SAPK2b2	Kinase	E. coli	GST	Human	71 kDa	14-244
p38 $\gamma$ /SAPK3, active	p38 $\gamma$ /SAPK3	Kinase	E. coli	GST	Human	69 kDa	14-246
p38 $\delta$ /SAPK4, active	p38 $\delta$ /SAPK4	Kinase	E. coli	GST	Human	68.9 kDa	14-249
p53 (expressed in E.coli)	p53	Kinase	E. coli	GST, c-Myc	Human	72 kDa	14-952
p70 S6 Kinase, active	p70S6K	Kinase	Sf21 insect cells	His6	Human	48.7 kDa	14-486
PAK1, active	PAK1	Kinase	Sf21 insect cells	FLAG	Human	46 kDa	14-927
PAK2, active	PAK2	Kinase	E. coli	His6	Human	62.8 kDa	14-481
PAK3, active	PAK3	Kinase	Sf21 insect cells	His6	Human	63.2 kDa	14-683
PAK4, active	PAK4	Kinase	Sf21 insect cells	GST, His6	Human	61.6 kDa	14-584
PAK5, active	PAK5	Kinase	Sf21 insect cells	His6	Human	37.4 kDa	14-699
PAK6, active	PAK6	Kinase	Sf21 insect cells	His6	Human	37.9 kDa	14-633
PAR-1B $\alpha$ , active	PAR-1B $\alpha$	Kinase	Sf21 insect cells	His6	Human	81.6 kDa	14-544
PASK, active	PASK	Kinase	Sf21 insect cells	His6	Human	41.5 kDa	14-701
PDGFR $\alpha$ (S50-end, V561D), active	PDGFR $\alpha$	Kinase	Sf21 insect cells	His6	Human	63.5 kDa	14-735
PDGFR $\alpha$ (S50-end,D842V), active	PDGFR $\alpha$	Kinase	Sf21 insect cells	His6	Human	63.5 kDa	14-729
PDGFR $\alpha$ , active	PDGFR $\alpha$	Kinase	Sf21 insect cells	His6	Human	63.5 kDa	14-467
PDGFR $\beta$ , active	PDGFR $\beta$	Kinase	Sf21 insect cells	His6	Human	63.7 kDa	14-463
PDHK2, active	PDHK2	Kinase	Sf21 insect cells	GST	Human	73 kDa	16-038-K
PDHK4, active	PDHK4	Kinase	Sf21 insect cells	GST	Human	74 kDa	15-024
PDK1, active	PDK1	Kinase	Sf21 insect cells	His6	Human	59 kDa	14-452
PEK, active	PEK	Kinase	Sf21 insect cells	His6	Human	71 kDa	14-916
PhK $\gamma$ 1, active	PhK $\gamma$ 1	Kinase	Sf21 insect cells	GST	Human	72 kDa	16-018-K
PhK $\gamma$ 2, active	PhK $\gamma$ 2	Kinase	E. coli	GST	Human	61.37 kDa	14-698
PI3 Kinase (p110 $\alpha$ (H1047R)/p85 $\alpha$ ) mouse	PI3K $\alpha$ /p85 $\alpha$	Kinase	Sf21 insect cells	His6	Mouse	129 kDa (p110 $\alpha$ H1047R), 83.6 kDa (p85 $\alpha$ )	14-787
PI3 Kinase (p110 $\alpha$ (E542K)/p85 $\alpha$ )	PI3K $\alpha$ /p85 $\alpha$	Kinase	Sf21 insect cells	His6	Human	125.3 kDa (p110 $\alpha$ E542K), 83.7 kDa (p85 $\alpha$ )	14-782
PI3 Kinase (p110 $\alpha$ (E545K)/p85 $\alpha$ )	PI3K $\alpha$ /p85 $\alpha$	Kinase	Sf21 insect cells	His6	Human	125.3 kDa (p110 $\alpha$ E542K), 83.7 kDa (p85 $\alpha$ )	14-783
PI3 Kinase (p110 $\alpha$ (E545K)/p85 $\alpha$ ), mouse	PI3K $\alpha$ /p85 $\alpha$	Kinase	Sf21 insect cells	His6	Mouse	129 kDa (p110 $\alpha$ E545K), 83.6 kDa (p85 $\alpha$ )	14-781
PI3 Kinase (p110 $\alpha$ /p65 $\alpha$ ), mouse	PI3K $\alpha$ /p65 $\alpha$	Kinase	Sf21 insect cells	His6	Mouse	129 kDa (p110 $\alpha$ ), 65.9 kDa (p65 $\alpha$ )	14-786
PI3 Kinase (p110 $\alpha$ /p65 $\alpha$ ), mouse	PI3K $\alpha$ /p65 $\alpha$	Kinase	Sf21 insect cells	His6	Mouse	125.3 kDa (p110 $\alpha$ ), 66 kDa (p65 $\alpha$ )	14-790
PI3 Kinase (p110 $\alpha$ /p85 $\alpha$ ) mouse	PI3K $\alpha$ /p85 $\alpha$	Kinase	Sf21 insect cells	His6	Mouse	129 kDa (p110 $\alpha$ ), 83.6 kDa (p85 $\alpha$ )	14-785

Recombinant Kinases and Phosphatases		Multiple sizes including bulk sizes are available					
Product Name	Target	Target Type	Expression System	Tag	Species	Molecular Weight	Base Catalog No.
PI3 Kinase (p110α/p85α), active	PI3Kα/p85α	Kinase	Sf21 insect cells	His6	Human	125.3 kDa (p110α), 83.7 kDa (p85α)	14-602
PI3 Kinase (p110β/p85α), active	PI3Kβ/p85α	Kinase	Sf21 insect cells	His6	Human	124 kDa (p110β), 83.7 kDa (p85α)	14-603
PI3 Kinase (p110β/p85β) mouse	PI3Kβ/p85β	Kinase	Sf21 insect cells	His6	Mouse	125.5 kDa (p110β), 81.4 kDa (p85β)	14-788
PI3 Kinase (p110δ/p85α), active	PI3Kδ/p85α	Kinase	Sf21 insect cells	His6	Human	121.6 kDa (p110δ), 83.7 kDa (p85α)	14-604
PI3 Kinase (p120γ)	PI3Kγ	Kinase	Sf21 insect cells	His6	Human	130 kDa	14-558
PI3K-C2α	PI3K-C2α	Kinase	Sf21 insect cells	His6	Human	161 kDa	14-906
PI3K-C2β	PI3K-C2β	Kinase	Sf21 insect cells	His6	Human	189 kDa	14-907
PI3K-C2γ	PI3K-C2γ	Kinase	Sf21 insect cells	GST	Human	112 kDa	14-910
PI3-Kinase (p110α(E542K)/p85α) mouse	PI3Kα/p85α	Kinase	Sf21 insect cells	His6	Mouse	129 kDa (p110α E542K), 84 kDa (p85α)	14-791
PI3-Kinase (p110α(H1047R)/p85α)	PI3Kα/p85α	Kinase	Sf21 insect cells	His6	Human	125 kDa (p110α H1047R), 84 kDa (p85α)	14-792
PI3-Kinase (p110β/p85α) mouse	PI3Kβ/p85α	Kinase	Sf21 insect cells	His6	Mouse	126 kDa (p110β), 84 kDa (p85α)	14-794
PI3-Kinase (p110δ/p85α) mouse	PI3Kδ/p85α	Kinase	Sf21 insect cells	His6	Mouse	124 kDa (p110δ), 84 kDa (p85α)	14-789
PI4KIIIa	PI4KIIIa	Kinase	Sf21 insect cells	FLAG	Human	233 kDa	14-908
PIK3C3, active	PIK3C3	Kinase	Sf21 insect cells	His6	Human	105 kDa	14-940
Pim1, active	Pim-1	Kinase	E. coli	GST	Human	62 kDa	14-573
PIM2, active	Pim-2	Kinase	E. coli	GST	Human	63.4 kDa	14-607
Pim-3, active	Pim-3	Kinase	Sf21 insect cells	His6	Human	39.6 kDa	14-738
PIP4K2α, active	PIP4K2α	Kinase	Sf21 insect cells	His6	Human	51 kDa	14-901
PIP5K1a, active	PIP5K1a	Kinase	Sf21 insect cells	His6	Human	60 kDa	14-844
PIP5K1γ, active	PIP5K1γ	Kinase	Sf21 insect cells	His6	Human	77 kDa	14-845
PKA, catalytic subunit, recombinant	PKA	Kinase	E. coli		Human	40.7 kDa	14-440
PKAcβ, active	PKAcβ	Kinase	Sf21 insect cells	His6	Human	45 kDa	15-007-K
PKC iota, active	PKCι	Kinase	Sf21 insect cells	His6	Human	68.6 kDa	14-505
PKC ζ, active	PKCζ	Kinase	Sf21 insect cells	His6	Human	68.9 kDa	14-525
PKC η, active	PKCη	Kinase	Sf21 insect cells	His6	Human	78.6 kDa	14-497
PKC θ, active	PKCθ	Kinase	Sf21 insect cells	His6	Human	83 kDa	14-444
PKCα, active	PKCα	Kinase	Sf21 insect cells	His6	Human	78 kDa	14-484
PKCβI, active	PKCβI	Kinase	Sf21 insect cells	His6	Human	78 kDa	14-503
PKCβII, active	PKCβII	Kinase	Sf21 insect cells	His6	Human	78.2 kDa	14-496
PKCγ, active	PKCγ	Kinase	Sf21 insect cells	His6	Human	79.5 kDa	14-483
PKCδ, active	PKCδ	Kinase	Sf21 insect cells	His6	Human	78.7 kDa	14-504
PKCε, active	PKCε	Kinase	Sf21 insect cells	His6	Human	84.6 kDa	14-518
PKCμ, active	PKCμ	Kinase	Sf21 insect cells	His6	Human	105 kDa	14-508
PKD2, active	PKD2	Kinase	Sf21 insect cells	His6	Human	100 kDa	14-506
PKD3, active	PKD3	Kinase	Sf21 insect cells	GST	Human	128 kDa	15-013-K
PKG1α, active	PKG1α	Kinase	Sf21 insect cells	His6	Human	80 kDa	14-688
PKG1β, active	PKG1β	Kinase	Sf21 insect cells	His6	Human	81.6 kDa	14-650
PKR, active	PKR	Kinase	Sf21 insect cells	GST, 10His	Human	64 kDa	14-955
Plk1, active	Plk1	Kinase	Sf21 insect cells	His6	Human	69.3 kDa	14-777
PLK3, active	PLK3	Kinase	Sf21 insect cells	His6	Human	36 kDa	14-572
Plk4, active	Plk4	Kinase	Sf21 insect cells	His6	Human	35 kDa	16-026-K
PP1α	PP1α	Phosphatase	E. coli		Human	37.6 kDa	14-595
PP5, active	PP5	Phosphatase	E. coli	GST	Human	83.8 kDa	14-778
PRAK, active	PRAK	Kinase	Sf21 insect cells	His6	Human	55 kDa	14-334
PRAK, unactive	PRAK	Kinase	Sf21 insect cells	His6	Human	54 kDa	14-335
PRK1, active	PRK1	Kinase	Sf21 insect cells	GST	Human	77 kDa	15-029-K
PRK2, active	PRK2	Kinase	Sf21 insect cells	His6	Human	59.9 kDa	14-549
PRKG2, active	PRKG2	Kinase	Sf21 insect cells	GST	Human	115 kDa	15-006-K
PRKX, active	PrKX	Kinase	Sf21 insect cells	His6	Human	44.7 kDa	14-677
PRP4, active	PRP4	Kinase	Sf21 insect cells	His6	Human	44 kDa	16-005-K
PTK5, active	PTK5	Kinase	Sf21 insect cells	His6	Human	38 kDa	14-693
PTP-1B	PTP-1B	Phosphatase	E. coli	GST	Human	64.1 kDa	14-621
PTPMEG-1	PTPMEG-1	Phosphatase	E. coli	GST	Human	83.1 kDa	14-642
PTPMEG-2	PTPMEG-2	Phosphatase	E. coli	GST	Human	62.5 kDa	14-592
PTPN22, active	PTPN22	Phosphatase	E. coli	GST	Human	63.3 kDa	14-768
PTPβ, active	PTPβ	Phosphatase	E. coli	GST	Human	69 kDa	14-948
Pyk2, active	Pyk2	Kinase	Sf21 insect cells	His6	Human	117.2 kDa	14-567
Raf-1 (truncated), active	c-RAF	Kinase	Sf21 insect cells	GST	Human	65 kDa	14-352
Ret (V804L), active	Ret	Kinase	Sf21 insect cells	GST	Human	79.2 kDa	14-758
Ret (V804M), active	Ret	Kinase	Sf21 insect cells	GST	Human	79.2 kDa	14-760
Ret, active	Ret	Kinase	Sf21 insect cells	GST	Human	79.2 kDa	14-570
RIPK1, active	RIPK1	Kinase	Sf21 insect cells	His6	Human	37 kDa	16-022-K
RIPK2, active	RIPK2	Kinase	Sf21 insect cells	His6	Human	37.8 kDa	14-612
ROKα/ROCK-II, active	ROKα/ROCK-II	Kinase	Sf21 insect cells	His6	Human	63.3 kDa	14-451
ROKα/ROCK-II, active, rat	ROKα/ROCK-II	Kinase	Sf21 insect cells	His6	Rat	66 kDa	14-338
ROKβ/ROCK-I, active	ROKβ/ROCK-I	Kinase	Sf21 insect cells	His6	Human	61.4 kDa	14-601
Ron, active	Ron	Kinase	Sf21 insect cells	GST	Human	74.2 kDa	14-581
Ros, active	Ros	Kinase	Sf21 insect cells	His6	Human	53.4 kDa	14-527
RPTPμ, active	RPTPμ	Phosphatase	E. coli	GST	Human	61.9 kDa	14-780
Rse, active	Rse	Kinase	Sf21 insect cells	His6	Human	50.4 kDa	14-535
Rsk1/MAPKAP Kinase 1a, active	Rsk1	Kinase	Sf21 insect cells	His6	Human	84 kDa	14-509
Rsk1/MAPKAP Kinase 1a, active, rat	Rsk1	Kinase	Sf21 insect cells	His6	Rat	88 kDa	14-479
Rsk2/MAPKAP Kinase 1b, active	Rsk2	Kinase	Sf21 insect cells	His6	Human	88 kDa	14-480
Rsk3, active	Rsk3	Kinase	Sf21 insect cells	His6	Human	87 kDa	14-462
Rsk4, active	Rsk4	Kinase	Sf21 insect cells	His6	Human	85.2 kDa	14-702
SAPK4, unactive	SAPK4	Kinase	E. coli	GST	Human	69 kDa	14-250
SBK1, active	SBK1	Kinase	Sf21 insect cells	GST	Human	74 kDa	16-029-K
SGK1 (S422D), unactive	SGK1	Kinase	Sf21 insect cells	His6	Human	48 kDa	14-332
SGK1 (61-59, S422D), active	SGK1	Kinase	Sf21 insect cells	His6	Human	48 kDa	14-331
SGK2 (hu, S4-end, S416D), unactive	SGK2	Kinase	Sf21 insect cells	His6	Human	45.7 kDa	14-636
SGK3, active	SGK3	Kinase	Sf21 insect cells	His6	Human	46.7 kDa	14-647
SGK3, unactive	SGK3	Kinase	Sf21 insect cells	His6	Human	46.8 kDa	14-648
SHP-1	SHP-1	Phosphatase	E. coli	GST	Human	95 kDa	14-591

Recombinant Kinases and Phosphatases					Multiple sizes including bulk sizes are available			
Product Name	Target	Target Type	Expression System	Tag	Species	Molecular Weight	Base Catalog No.	
SHP-2	SHP-2	Phosphatase	E. coli	GST	Human	63.7 kDa	14-622	
SIK, active	SIK	Kinase	Sf21 insect cells	His6	Human	36.2 kDa	14-652	
SIK2, active	SIK2	Kinase	Sf21 insect cells	GST	Human	59 kDa	15-010	
SIK3, active	SIK3	Kinase	Sf21 insect cells	GST, His6	Human	63 kDa	15-011	
SLK, active	SLK	Kinase	Sf21 insect cells	His6	Human	46 kDa	15-027-K	
SNRK, active	SNRK	Kinase	Sf21 insect cells	His6	Human	58 kDa	14-932	
Src (1-530), active	Src	Kinase	Sf21 insect cells	His6	Human	61.7 kDa	14-746	
Src (T341M), active	Src	Kinase	Sf21 insect cells	His6	Human	61.7 kDa	14-748	
Src, active	Src	Kinase	Sf21 insect cells	His6	Human	61.7 kDa	14-326	
SRMS, active	SRMS	Kinase	Sf21 insect cells	GST	Human	58 kDa	16-036-K	
SRPK1, active	SRPK1	Kinase	E. coli	GST	Human	100 kDa	14-564	
SRPK2, active	SRPK2	Kinase	E. coli	GST	Human	100 kDa	14-666	
STK16, active	STK16	Kinase	Sf21 insect cells	GST, His6	Human	62 kDa	16-035-K'	
STK25, active	STK25	Kinase	Sf21 insect cells	GST, His6	Human	63 kDa	14-929	
STK32A, active	STK32A	Kinase	Sf21 insect cells	His6	Human	50 kDa	16-011	
STK32B, active	STK32B	Kinase	Sf21 insect cells	His6	Human	53 kDa	16-021-K	
STK32C, active	STK32C	Kinase	Sf21 insect cells	His6	Human	59 kDa	16-012-K	
STK33, active	STK33	Kinase	Sf21 insect cells	His6	Human	62.4 kDa	14-671	
STK39, active	STK39	Kinase	Sf21 insect cells	GST	Human	91 kDa (STK39), 67 kDa (MO25α)	16-049	
Syk, active	Syk	Kinase	Sf21 insect cells	His6	Human	73 kDa	14-314	
TAF1L, active	TAF1L	Kinase	Sf21 insect cells	His6	Human	49 kDa	16-024-K	
TAK1-TAB1 fusion, active	Tak1	Kinase	Sf21 insect cells	His6	Human	42.8 kDa	14-600	
TAO1, active	TAO1	Kinase	Sf21 insect cells	His6	Human	41 kDa	14-749	
TAO2, active	TAO2	Kinase	Sf21 insect cells	His6	Human	138.3 kDa	14-736	
TAO3, active	TAO3	Kinase	Sf21 insect cells	His6	Human	50.2 kDa	14-745	
TBK1, active	TBK1	Kinase	Sf21 insect cells	His6	Human	85.7 kDa	14-628	
TCPTP	TCPTP	Phosphatase	E. coli		Human	39.7 kDa	14-646	
Tec, activated	Tec	Kinase	Sf21 insect cells	His6	Human	57 kDa	14-801	
TGFBR-1, active	TGFBR1	Kinase	Sf21 insect cells	GST	Human	62 kDa	14-912	
TGFBR2, active	TGFBR2	Kinase	Sf21 insect cells	His6	Human	45 kDa	16-014-K	
Tie2 (R849W), active	Tie2	Kinase	Sf21 insect cells	His6	Human	42 kDa	14-763	
Tie2 (Y1108F), active	Tie2	Kinase	Sf21 insect cells	His6	Human	42 kDa	14-766	
Tie2 (Y897S), active	Tie2	Kinase	Sf21 insect cells	His6	Human	42 kDa	14-764	
Tie2, active	Tie2	Kinase	Sf21 insect cells	His6	Human	42 kDa	14-540	
TLK1, active	TLK1	Kinase	Sf21 insect cells	His6	Human	91 kDa	14-938	
TLK2, active	TLK2	Kinase	Sf21 insect cells	His6	Human	48.8 kDa	14-739	
TMDP	TMDP	Phosphatase	E. coli	GST	Human	49 kDa	14-767	
TNIK, active	TNIK	Kinase	Sf21 insect cells	GST	Human	69 kDa	15-016-K	
TRB2, active	TRB2	Kinase	Sf21 insect cells	His6	Human	43 kDa	16-016-K	
TrkA, active	TrkA	Kinase	Sf21 insect cells	His6	Human	41 kDa	14-571	
TrkB, active	TrkB	Kinase	Sf21 insect cells	His6	Human	42.9 kDa	14-507	
TrkC, active	TrkC	Kinase	Sf21 insect cells	GST, His6	Human	64 kDa	14-922	
TSSK1, active	TSSK1	Kinase	Sf21 insect cells	His6	Human	46 kDa	14-670	
TSSK2, active	TSSK2	Kinase	Sf21 insect cells	His6	Human	44.8 kDa	14-632	
TSSK3, active	TSSK3	Kinase	Sf21 insect cells	His6	Human	31 kDa	15-021	
TSSK4, active	TSSK4	Kinase	Sf21 insect cells	GST	Human	65 kDa	15-017	
TTBK1, active	TTBK1	Kinase	Sf21 insect cells	GST, His6	Human	82 kDa	15-018-K	
TTBK2, active	TTBK2	Kinase	Sf21 insect cells	GST	Human	65 kDa	15-019	
TTK, active	TTK	Kinase	Sf21 insect cells	GST	Human	124 kDa	15-014	
Txk, active	Txk	Kinase	Sf21 insect cells	His6	Human	35.3 kDa	14-761	
TYK2, active	TYK2	Kinase	Sf21 insect cells	His6	Human	40 kDa	14-924	
ULK1, active	ULK1	Kinase	Sf21 insect cells	His6	Human	39 kDa	14-959	
ULK2, active	ULK2	Kinase	Sf21 insect cells	GST, His6	Human	62 kDa	14-772	
ULK3, active	ULK3	Kinase	Sf21 insect cells	His6	Human	62 kDa	14-755	
VHR	VHR	Phosphatase	E. coli	GST	Human	47.8 kDa	14-594	
VRK1, active	VRK1	Kinase	Sf21 insect cells	His6	Human	50 kDa	16-033-K	
VRK2, active	VRK2	Kinase	Sf21 insect cells	His6	Human	43 kDa	14-732	
Wee1, active	WEE1	Kinase	Sf21 insect cells	GST, His6	Human	77 kDa	14-925	
Wee1B, active	Wee1B	Kinase	Sf21 insect cells	His6	Human	67 kDa	16-006-K	
WNK1, active	WNK1	Kinase	Sf21 insect cells	His6	Human	58 kDa	16-007-K	
WNK2, active	WNK2	Kinase	Sf21 insect cells	His6	Human	41.2 kDa	14-678	
WNK3, active	WNK3	Kinase	Sf21 insect cells	His6	Human	53.1 kDa	14-658	
WNK4, active	WNK4	Kinase	Sf21 insect cells	His6	Human	54 kDa	16-047-K	
Yes, active	Yes	Kinase	Sf21 insect cells	His6	Human	63 kDa	14-478	
YopH, active, Yersinia	YopH	Phosphatase	E. coli	GST	Yersinia	77.4 kDa	14-590	
ZAK, active	ZAK	Kinase	Sf21 insect cells	GST	Human	119 kDa	15-015	
ZAP-70, active	ZAP-70	Kinase	Sf21 insect cells	His6	Human	70 kDa	14-404	
ZIPK, active	ZIPK	Kinase	E. coli	GST	Human	64.3 kDa	14-608	





 eurofins

DiscoverX