

HTRF[®] cellular phospho-protein assays



PATHWAY-READOUT

PHYSIOLOGICALLY RELEVANT RESULTS FOR FAST FLOWING RESEARCH

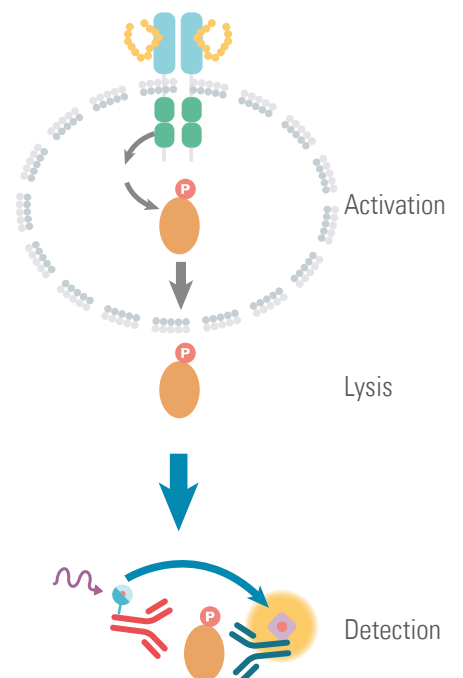
Time may be money, but relevant results are golden.

Cisbio's smart pathway readout tools simplify and accelerate your research projects, from basic biological investigations to high-throughput drug screening. They are ideal for both small molecule and biotherapeutic pharmacology.

Based on HTRF, Cisbio's unique Homogeneous Time-Resolved Fluorescence detection platform, this no-wash cell-based system offers add-and-read assays for physiologically relevant results, ideal for studying signaling pathways in the cell.

Progress faster, with more relevant cell signaling results using cell-based HTRF phospho-assays

- Greater specificity and high Z' scores for robust, reproducible results
- Fewer false negatives and repeat runs
- Adapts easily to all throughput requirements
- Simple wash free add-and-read protocol
- High species cross-reactivity



ORDERING INFORMATION

PRODUCT	PART#	PRODUCT	PART#	PRODUCT	PART#
4EBP1 phospho-T37/46 kit	64BP1PEG	GSK3 alpha total kit	64GTAPEG	RB phospho-S780 kit NEW	63ADK104PEG
Alpha Synuclein phospho-S129 kit	6FSYNPEG	GSK3 beta phospho-S9 kit	64GPBPEG	RB phospho-S807/811 kit NEW	63ADK105PEG
Alpha Synuclein total kit	6FNSYPEG	GSK3 beta total kit	64GTBPEG	RB total kit NEW	63ADK106PEG
ACC phospho-S79 kit	64ACCPPEG	H2AX phospho-S139 kit	64H2XPEG	S6RP phospho-S235/236 kit	64RP6PEG
Advanced ERK phospho-T202/Y204 kit	64AERPEG	HER2 phospho-Y1221/1222 kit	64HR2PEG	SHP1 phospho-Y564 kit	64SH1PEG
AKT1 phospho-S473 kit	64AKT078PEG	HER2 total kit	64NR2PEG	SHP1 total kit	64NH1PEG
AKT1 total kit	64AKT079PEG	HER3 phospho-Y1289 kit	64HR3PEG	SHP2 phospho-Y542 kit	64SH2PEG
AKT2 phospho-S473 kit	64AKT080PEG	HER3 total kit	64NR3PEG	SHP2 total kit	64NH2PEG
AKT2 total kit	64AKT081PEG	Histone H3 phospho-S10 kit	64HH3PEG	SLP-76 phospho-S376 kit	63ADK076PEG
AKT3 phospho-S473 kit	64AKT082PEG	Histone H3 phospho-T3 kit	63ADK061PEG	SLP-76 total kit	63ADK077PEG
AKT3 total kit	64AKT083PEG	Histone H3 total kit	64NH3PEG	SMAD1 phospho-S463/465 kit	63ADK062PEG
AKT phospho-S473 kit	64AKSPEG	IKKB phospho-S177/181 kit	64KKBPEG	SMAD1 total kit	63ADK063PEG
AKT phospho-T308 kit	64AKTPEG	IKKB total kit	63ADK097PEG	SMAD2 phospho-S465/467 kit	63ADK102PEG
AKT total kit	64NKTPEG	IRAK3 total kit	63ADK101PEG	SMAD2 total kit	63ADK103PEG
AMPK alpha phospho-T172 kit	64MPKPEG	IRF3 phospho-S385/386 kit	6FRF3PEG	SMAD3 phospho-S423/425 kit	63ADK025PEG
AMPK alpha total kit	63ADK060PEG	IR beta phospho-Y1150/1151 kit	63ADK016PEG	SMAD3 total kit	64ND3PEG
ATF2 phospho-T71 kit	63ADK015PEG	IR beta total kit	63ADK019PEG	c-Src phospho-Y419 kit	64SRCPEG
BAD phospho-S112 kit	64BADPEG	IRS1 phospho-S307(m)/S312(h) kit	64IRSPEG	c-Src total kit	64NSRPEG
Beta Catenin phospho-T41/S33/37 kit	64CATPEG	IRS1 total kit	64NRSPEG	STAT1 phospho-Y701 kit	63ADK026PEG
Beta Catenin total kit	64NCAPEG	IRS2 total kit	64IRTPEG	STAT1 total kit	63ADK096PEG
BTK phospho-Y223 kit	63ADK017PEG	JNK phospho-T183/Y185 kit	64JNKPEG	STAT3 phospho-Y705 kit	62AT3PEG
BTK total kit	63ADK064PEG	LRRK2 phospho-S935 kit	6FLRKPEG	STAT3 total kit	64NT3PEG
cJun phospho-S63 kit	64JUNPEG	LRRK2 total kit	6FNRKPEG	STAT5 phospho-Y694 kit	64AT5PEG
cRaf phospho-S338 kit	63ADK018PEG	MEK1 phospho-S218/222 kit	64ME1PEG	STAT6 phospho-S366 kit	64AT6PEG
CREB phospho-S133 kit	64CREPEG	MEK1 phospho-S298 kit	63ADK021PEG	STING phospho-Y641 kit	64STGPEG
CREB total kit	63ADK052PEG	MEK1 total kit	64NE1PEG	STING total kit	64NTGPEG
EGFR phospho-pan kit	64HR1PEG	MEK1/2 phospho-S218/222 kit	64ME2PEG	Tau phospho-S202/T205 kit	64TS2PEG
EGFR phospho-Y1068 kit	64EG1PEG	MNK-1 phospho-T197/202 kit	63ADK022PEG	Tau phospho-S404 kit	64TS4PEG
EGFR total kit	64NG1PEG	mTOR phospho-S2448 kit	64TORPEG	TBK1 phospho-S172 kit	64TBKPEG
EIF4E phospho-S209 kit	64EF4PEG	NDRG1 phospho-S330 kit	63ADK093PEG	TBK1 total kit	64NTBPEG
ERK phospho-T202/Y204 kit	64ERKPEG	NDRG1 phospho-T346 kit	63ADK092PEG	ULK1 phospho-S757 kit	64ULKPEG
EIF2 alpha phospho-S52 kit	64EF2PEG	NDRG1 total kit	63ADK094PEG	VASP phospho-S157 kit	63ADK066PEG
EIF2 alpha total kit	64NEFPEG	NFkB phospho-S536 kit	64NFBPEG	VASP phospho-S239 kit	63ADK065PEG
ERK total kit	64NRKPEG	NFkB total kit	64NFTPEG	VASP total kit	63ADK067PEG
FAK phospho-Y397 kit	64FAKPEG	Phospholamban phospho-T17 kit	63ADK074PEG	VEGFR2 Phospho-Y1175	63ADK041PEG
FAK total kit	64NAKPEG	Phospholamban total kit	63ADK075PEG	YAP phospho-S127 kit	64YAPPEG
FGFR2 phospho-Y653/654 kit	63ADK054PEG	PDGFR beta phospho-Y751 kit	64PDGPEG	YAP total kit	64YATPEG
FGFR2 total kit	63ADK057PEG	P38 phospho-T180/Y182 kit	64P38PEG	ZAP-70 phospho-Y319 kit	64ZAPPEG
FoxO1 phospho-S256 kit	64FOXPEG	P53 phospho-S15 kit	64P53PEG	ZAP-70 total kit	64ZATPEG
FoxO1 total	64NFOPEG	P53 total kit	64T53PEG		
GSK3 alpha phospho-S21 kit	64GPAPEG	p70S6K phospho-T389 kit	64S6KPEG		

500 tests each, bulk size available

PRODUCT	PART#
Alpha-Tubulin Housekeeping kit	64ATUBPEG

500 tests each, bulk size available



EXPERT SERVICES TO SATISFY YOUR EXACT NEEDS

If you have specific requirements for quantifying cellular phospho-proteins or other signaling proteins.

Call us or visit www.cisbio.com

FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567 China +86 21 5018 9880
Japan +81-(0)43-306-8712 Visit www.cisbio.com to find a list of our regional distributors

