

# HTRF<sup>®</sup> cellular phospho-protein assays



PATHWAY-READOUT

## PHYSIOLOGICALLY RELEVANT RESULTS FOR FAST FLOWING RESEARCH

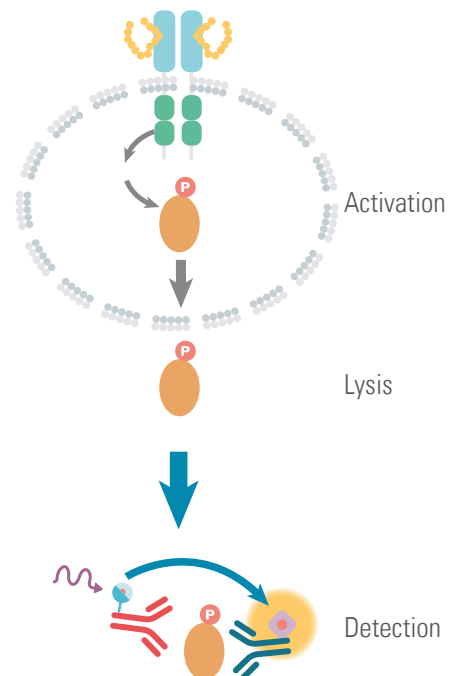
Time may be money, but relevant results are golden.

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	eNOS phospho-S1177 kit <b>NEW</b>	64ENOSS7PEG	MEK1 total kit	64NE1PEG
Alpha Synuclein phospho-S129 kit	6FSYNPEG	eNOS total kit <b>NEW</b>	64ENOSTPEG	MEK1/2 phospho-S218/222 kit	64ME2PEG
Alpha Synuclein total kit	6FNSTYPEG	ERK phospho-T202/Y204 kit	64ERKPEG	MKK4 phospho-S258 kit	64MK4PEG
ACC phospho-S79 kit	64ACCPEG	ERK Advanced phospho-T202/Y204 kit	64AERPEG	MKK4 total kit	64NK4PEG
ACC total kit	64NACPEG	ERK total kit	64NRKPEG	MNK-1 phospho-T197/202 kit	63ADK022PEG
AKT1 phospho-S473 kit	64AKT078PEG	FAK phospho-Y397 kit	64FAKPEG	mTOR phospho-S2448 kit	64TORPEG
AKT1 total kit	64AKT079PEG	FAK total kit	64NAKPEG	NDRG1 phospho-S330 kit	63ADK093PEG
AKT2 phospho-S473 kit	64AKT080PEG	FGFR2 phospho-Y653/654 kit	63ADK054PEG	NDRG1 phospho-T346 kit	63ADK092PEG
AKT2 total kit	64AKT081PEG	FGFR2 total kit	63ADK057PEG	NDRG1 total kit	63ADK094PEG
AKT3 phospho-S473 kit	64AKT082PEG	FoxO1 phospho-S256 kit	64FOXPEG	NFkB phospho-S536 kit	64NFBPEG
AKT3 total kit	64AKT083PEG	FoxO1 total kit	64NFOPEG	NFkB total kit	64NFTPEG
AKT phospho-S473 kit	64AKSPEG	GSK3 alpha phospho-S21 kit	64GPAPEG	Phospholamban phospho-S16 kit	64PLN16PEG
AKT phospho-T308 kit	64AKTPEG	GSK3 alpha total kit	64GTAPEG	Phospholamban phospho-T17 kit	63ADK074PEG
AKT total kit	64NKTPEG	GSK3 beta phospho-S9 kit	64GPBPEG	Phospholamban total kit	63ADK075PEG
AMPK alpha phospho-T172 kit	64MPKPEG	GSK3 beta total kit	64GTBPEG	PDGFR beta phospho-Y751 kit	64PDGPEG
AMPK alpha total kit	63ADK060PEG	H2AX phospho-S139 kit	64H2XPEG	PLCg1 phospho-Y783 kit <b>NEW</b>	64PLCG1Y3PEG
ATF2 phospho-T71 kit	63ADK015PEG	HER2 phospho-Y1221/1222 kit	64HR2PEG	PLCg1 total kit <b>NEW</b>	64PLCG1TPEG
BAD phospho-S112 kit	64BADPEG	HER2 total kit	64NR2PEG	PLCg2 phospho-Y1214 kit	64PLCG2Y7PEG
Beta Catenin phospho-T41/S33/37 kit	64CATPEG	HER3 phospho-Y1289 kit	64HR3PEG	PLCg2 total kit	64PLCG2TPEG
Beta Catenin total kit	64NCAPEG	HER3 total kit	64NR3PEG	P38 phospho-T180/Y182 kit	64P38PEG
BTK phospho-Y223 kit	63ADK017PEG	Histone H3 phospho-S10 kit	64HH3PEG	P53 phospho-S15 kit	64P53PEG
BTK total kit	63ADK064PEG	Histone H3 phospho-T3 kit	63ADK061PEG	P53 total kit	64T53PEG
CDK2 phospho-Y15 kit	64CDK2Y15PEG	Histone H3 total kit	64NH3PEG	p70S6K phospho-T389 kit	64S6KPEG
CDK2 total kit	64CDK2TPEG	IKKB phospho-S177/181 kit	64KKBPEG	RB phospho-S780 kit	63ADK104PEG
CDK4 total kit	64CDK4TPEG	IKKB total kit	63ADK097PEG	RB phospho-S807/811 kit	63ADK105PEG
CDK6 total kit	64CDK6TPEG	IRAK3 total kit	63ADK101PEG	RB total kit	63ADK106PEG
Cereblon total kit <b>NEW</b>	64CRBNPEG	IRF3 phospho-S385/386 kit	6FRF3PEG	S6RP phospho-S235/236 kit	64RP6PEG
cJun phospho-S63 kit	64JUNPEG	IR beta phospho-Y1150/1151 kit	63ADK016PEG	SHP1 phospho-Y564 kit	64SH1PEG
cRaf phospho-S338 kit	63ADK018PEG	IR beta total kit	63ADK019PEG	SHP1 total kit	64NH1PEG
CREB phospho-S133 kit	64CREPEG	IRS1 phospho-S307(m)/S312(h) kit	64IRSPEG	SHP2 phospho-Y542 kit	64SH2PEG
CREB total kit	63ADK052PEG	IRS1 total kit	64NRSPEG	SHP2 total kit	64NH2PEG
EGFR phospho-pan kit	64HR1PEG	IRS2 total kit	64IRTPEG	SLP-76 phospho-S376 kit	63ADK076PEG
EGFR phospho-Y1068 kit	64EG1PEG	JNK phospho-T183/Y185 kit	64JNKPEG	SLP-76 total kit	63ADK077PEG
EGFR total kit	64NG1PEG	LRRK2 phospho-S935 kit	6FLRKPEG	SLP-76 mouse phospho-S376 kit	63ADK109PEG
EIF4E phospho-S209 kit	64EF4PEG	LRRK2 total kit	6FNKPEG	SLP-76 mouse total kit	63ADK110PEG
EIF2 alpha phospho-S52 kit	64EF2PEG	MEK1 phospho-S218/222 kit	64ME1PEG	SMAD1 phospho-S463/465 kit	63ADK062PEG
EIF2 alpha total kit	64NEFPEG	MEK1 phospho-S298 kit	63ADK021PEG	SMAD1 total kit	63ADK063PEG

PRODUCT	PART#	PRODUCT	PART#	PRODUCT	PART#
SMAD2 phospho-S465/467 kit	63ADK102PEG	SYK phospho-Y525 kit	64SYKY525PEG	VEGFR2 Phospho-Y1175	63ADK041PEG
SMAD2 total kit	63ADK103PEG	SYK total kit	64SYKTPEG	YAP phospho-S127 kit	64YAPPEG
SMAD3 phospho-S423/425 kit	63ADK025PEG	Tau phospho-S202/T205 kit	64TS2PEG	YAP total kit	64YATPEG
SMAD3 total kit	64ND3PEG	Tau phospho-356 kit	64TS356PEG	ZAP-70 phospho-Y319 kit	64ZAPPEG
c-Src phospho-Y419 kit	64SRCPEG	Tau phospho-S404 kit	64TS4PEG	ZAP-70 total kit	64ZATPEG
c-Src total kit	64NSRPEG	Tau phospho-S422 kit	64TS422PEG		
STAT1 phospho-Y701 kit	63ADK026PEG	Tau phospho-T181 kit	64TT181PEG		
STAT1 total kit	63ADK096PEG	Tau total kit	64NTAUPEG		
STAT3 phospho-Y705 kit	62AT3PEG	TBK1 phospho-S172 kit	64TBKPEG		
STAT3 total kit	64NT3PEG	TBK1 total kit	64NTBPEG		
STAT5 phospho-Y694 kit	64AT5PEG	ULK1 phospho-S757 kit	64ULKPEG		
STAT6 phospho-S366 kit	64AT6PEG	VASP phospho-S157 kit	63ADK066PEG		
STING phospho-Y641 kit	64STGPEG	VASP phospho-S239 kit	63ADK065PEG		
STING total kit	64NTGPEG	VASP total kit	63ADK067PEG		

*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](http://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 10 8080 9288 Japan +81-(0)43-306-8712

Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors



cisbio