

HTRF[®] cell based phospho-assays

1 x 96 tests



PATHWAY-READOUT

HAVE YOU EVER WANTED TO RUN 96 WESTERN BLOTS IN 1 DAY?

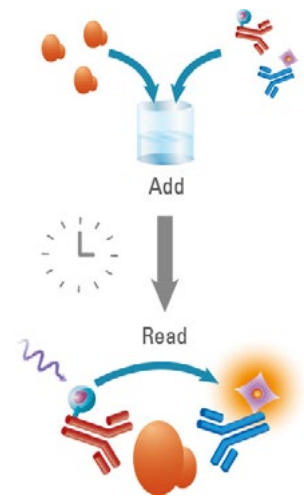
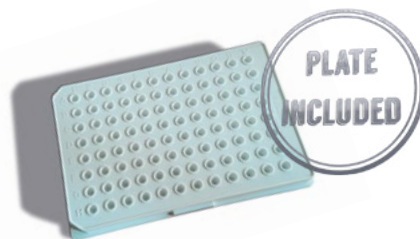
Discover Cisbio's straightforward protocols for cell-based phospho protein assays:

- 16 μ L of sample (cell lysate, tissue homogenate ...)
- 4 μ L of detection reagents

Just add, incubate and read!

These kits contain everything you need to get started:

- Detection antibodies, buffers, control, protocol
- 96-well low volume plate



HTRF[®] technology, an effective alternative to ELISA & Western Blot

- Simple wash-free add-and-read protocol saves you time
- Low consumption saves on precious samples

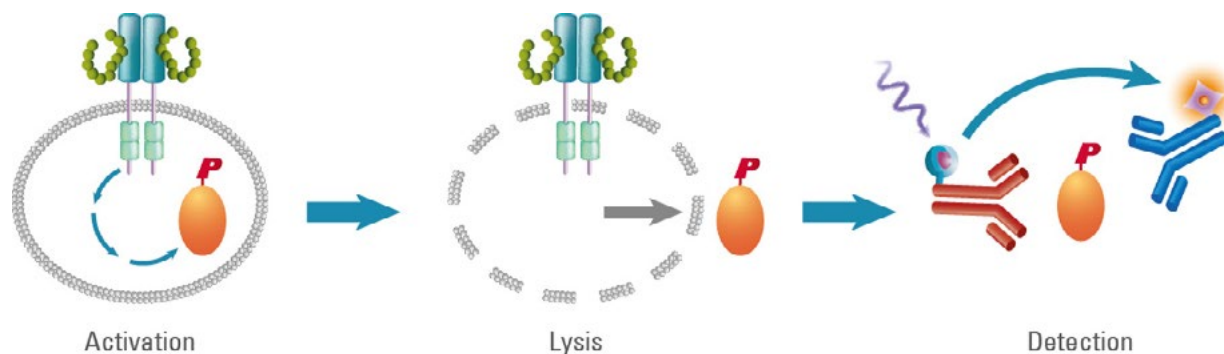
Check out a protocol video: www.cisbio.com/ERK-video

cisbio
INTERACTION IS EVERYTHING

HOW IT WORKS

AN ASSAY PROTOCOL THAT IS SIMPLER AND MORE FLEXIBLE

After lysis of the cell membrane, phosphorylated proteins and total protein populations can be detected using HTRF kit reagents and most TR-FRET multimode plate readers.



PART NUMBER#	PRODUCT DESCRIPTION	TESTS
64ACCPET	ACC phospho-S79 kit	96 tests
64AKSPET	AKT phospho-S473 kit	96 tests
64NKTPET	AKT total kit	96 tests
64MPKPET	AMPK phospho-T172 kit	96 tests
64BADPET	BAD phospho-S112 kit	96 tests
64CREPET	CREB phospho-S133 kit	96 tests
64AERPET	Advanced ERK phospho-T202 /Y204 kit	96 tests
64NRKPET	ERK total kit	96 tests
64NFBPET	NFkB phospho-S536 kit	96 tests
64P38PET	P38 phospho-T180/Y182 kit	96 tests
64P53PET	P53 phospho-S15 kit	96 tests
64T53PET	P53 total kit	96 tests
63ADK076PET	SLP-76 phospho-S376 kit	96 tests
63ADK062PET	SMAD1 phospho-S463/465 kit	96 tests
62AT3PET	STAT3 phospho-Y705 kit	96 tests

This list is updated regularly with new references.

EXPERT SERVICES TO SATISFY YOUR PRECISE NEEDS



In addition to providing ready-to-use kits with all the necessary reagents, Cisbio's team of experienced experts works hand-in-hand with our clients to create custom solutions. When you have specific requirements for quantifying cellular phospho-proteins or other signaling proteins, Cisbio is the partner you can trust to help develop the assays you need.

Keep your therapeutic research moving forward with faster, more robust, physiologically relevant results. Call us or visit www.cisbio.com/cellular-kinase-assays

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 Cisbio.com to find a list of our regional distributors

