

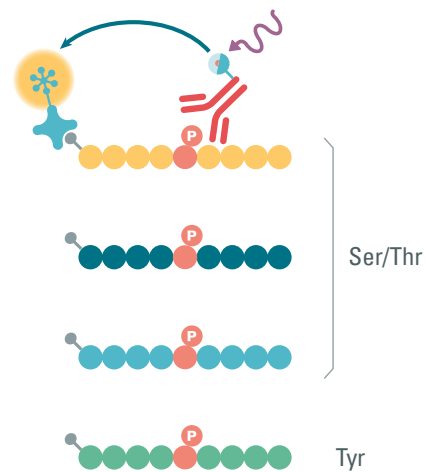
HTRF® KINEASE™ TYR AND SER/THR KITS



A UNIVERSAL ASSAY PLATFORM FOR SCREENING SERINE/THREONINE AND TYROSINE KINASE ACTIVITY

Serine/Threonine & Tyrosine kinases are major targets in drug discovery. Cisbio has combined its HTRF technology and its expertise in the kinase field to design a universal and sensitive assay packed into ready-to-use kits for measuring the activity of large panels of Serine/Threonine and Tyrosine kinases. Simple and fast—ideal for profiling and HTS. Over 200 different kinases have been validated using HTRF KinEASE kits.

- Low enzyme consumption
- No limitation in ATP concentration
- Results in under 2 hours
- Easy miniaturization to <4 µL
- Discovery kit for Ser/Thr kinases
- One universal substrate for Tyr kinases



Four KinEASE substrates cover a broad selection of Ser/Thr and Tyr kinases.

KINASES

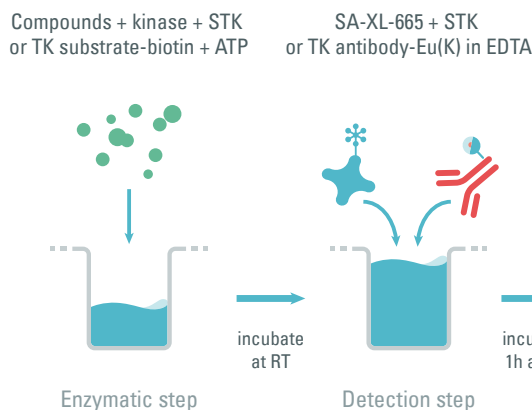
ASSAY PROTOCOL

HTRF KinEASE assays use a biotinylated substrate, a phospho-specific monoclonal antibody labeled with Eu³⁺ cryptate (donor) and XL665 labeled streptavidin (acceptor). After phosphorylation, the anti-phospho monoclonal antibody will recognize the site with its phosphorylation mark and the proximity of donor and acceptor will consequently lead to a TR-FRET signal.

The basic assay involves two steps: the enzymatic step followed by detection.

Enzymatic step: The kinase is incubated in the presence or absence of compounds and the appropriate substrate. ATP is added to start the reaction.

Detection step: Eu³⁺ cryptate and XL665 conjugate are prepared in the supplemented detection buffer and added to the reaction. The EDTA in the detection buffer will stop the enzymatic reaction. The two detection reagents can be pre-mixed and added in a single dispensing step providing enhanced speed and simplicity.



ORDERING INFORMATION:

PRODUCT	TESTS	PART #	COMPANION PRODUCTS	PART #
HTRF KinEASE-STK Discovery kit*	1,000	62ST0PEB	SEB-Suppl. Enzymatic Buffer—12.5 nmoles	61SEBALB
HTRF KinEASE-STK S1 kit	1,000	62ST1PEB	STK-Substrate 1-biotin—500 µg	61ST1BLC
HTRF KinEASE-STK S1 kit	20,000	62ST1PEC	STK-Substrate 1-biotin—50 µg	61ST1BLE
HTRF KinEASE-STK S1 kit	100,000	62ST1PEJ	STK-Substrate 2-biotin—500 µg	61ST2BLC
HTRF KinEASE-STK S2 kit	1,000	62ST2PEB	STK-Substrate 2-biotin—50 µg	61ST2BLE
HTRF KinEASE-STK S2 kit	20,000	62ST2PEC	STK-Substrate 3-biotin—500 µg	61ST3BLC
HTRF KinEASE-STK S2 kit	100,000	62ST2PEJ	STK-Substrate 3-biotin—50 µg	61ST3BLE
HTRF KinEASE-STK S3 kit	1,000	62ST3PEB	TK Substrate-biotin—500 µg	61TK0BLC
HTRF KinEASE-STK S3 kit	20,000	62ST3PEC	TK Substrate-biotin—50 µg	61TK0BLE
HTRF KinEASE-STK S3 kit	100,000	62ST3PEJ	Enzymatic KinEASE buffer 5X—10 mL	62EZBFDC
HTRF KinEASE-TK kit	1,000	62TK0PEB	Enzymatic KinEASE buffer 5X—50 mL	62EZBFDD
HTRF KinEASE-TK kit	20,000	62TK0PEC	HTRF KinEASE detection buffer—40 mL	62SDBRDD
HTRF KinEASE-TK kit	100,000	62TK0PEJ	HTRF KinEASE detection buffer—200 mL	62SDBRDF

*For screening the activity of Ser/Thr Kinases, Cisbio offers a special Discovery kit, which includes all three Ser/Thr substrates for you to easily determine and choose the optimal setup for your assay.

For more information, please visit www.htrf.com/kinases-cell-signaling

INTERACTION IS EVERYTHING

From the chemistry of our assays to the collaborative nature of our relationship with customers, Cisbio is founded on the belief that “interaction is everything”. Talk to us. And discover a partner dedicated to providing exactly what you need:

- Convenient ready-to-use kits
- Custom assay capabilities
- Expert technical support
- Complete confidentiality
- Trusted and proven HTRF technology

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 to find a list of our regional distributors



cisbio