



HTRF® KinEASE™

A universal expanded platform to address Serine/Threonine & Tyrosine kinases

APPLICATION NOTE

ABSTRACT Kinases are enzymes that play a central role in various signal transduction pathways involved in the control of cell growth, metabolism, differentiation and apoptosis. They exert their effect by catalysing the transfer of a phosphate group from ATP onto a target substrate (protein/peptide), which then becomes activated and performs a specific function.

HTRF KinEASE are universal tools for assessing Serine/Threonine (STK) and Tyrosine Kinase (TK) activity. A typical development for an HTRF KinEASE assay consists of five steps, and is described in this document.

Protein kinases are the second most important group of drug targets after GPCRs. The approval of two kinase-targeting drugs, Genentech's Herceptin® in 1998 and Novartis's Gleevec® in 2001, and their success in cancer therapeutics, confirm the trend towards screening new Protein Kinase Inhibitors (PKI) of various kinase targets for use in different therapeutic areas.

This growing interest in screening kinases has prompted the development of many assay technologies.

The new HTRF KinEASE TK kit for the measurement of Tyrosine Kinase (TK) activities combines a universal peptide substrate and a single proprietary monoclonal antibody with Cisbio's HTRF (Homogeneous Time-Resolved Fluorescence) technology, a highly sensitive and robust technology for the detection of molecular interactions of proteins in vitro.

HTRF KinEASE STK is the fourth kit in the HTRF KinEASE platform, developed in collaboration with Millipore (Upstate) for profiling and HTS of Serine/Threonine and Tyrosine Kinases. The platform's other three kits are dedicated to the Ser/Thr kinases group and combine three universal biotinylated substrates, S1, S2 and S3 and a monoclonal antibody. The new HTRF KinEASE kit contains a generic substrate and a monoclonal antibody.

HTRF KINEASE: ASSAY IMPLEMENTATION

HTRF KinEASE screenings limit assay development time and are easily miniaturizable and flexible, meaning the assay can be performed under a wide range of kinase assay conditions, for instance with low consumption of enzyme or with any ATP concentration. The assay is run in two main steps, the enzymatic (kinase reaction) step followed by the detection step with HTRF reagents (Fig. 1). The assay is started by the addition of ATP (Step 1) and stopped by the addition of the HTRF detection reagents containing EDTA (Step 2).

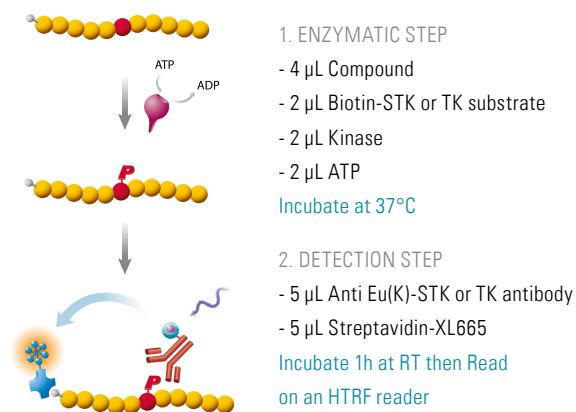


Fig. 1: HTRF KinEASE assay principle: All the assays based on HTRF KinEASE involve two steps, the enzymatic step and the detection step with HTRF reagents. Step 1: During the enzymatic step the substrate-biotin is incubated with the kinase of interest. ATP is added to start the reaction. Step 2: The detection reagent catches the phosphorylated substrate and the resulting TR-FRET is proportional to the phosphorylation level.

A STRAIGHTFORWARD KINASE ASSAY DEVELOPMENT

A typical development for an HTRF KinEASE assay consists of five steps, and is described in a document supplied with the kit. The conditions of each step are described for MAPKAP-K2 using HTRF KinEASE STK-S1.

1. ENZYME TITRATION

This step gives the optimal kinase concentration, i.e. that for which the signal reaches 80% of the maximum (EC_{80}). Kinase is used at concentrations ranging from 0.10 ng/well to 10 ng/well, and incubated 30 min with the substrate-biotin (1 μ M), and a non-limiting ATP concentration (100 μ M). The reaction is stopped by the addition of the HTRF detection reagents in EDTA. The substrate-biotin/SA-XL665 ratio of 8/1 (i.e. 62.2 nM SAXL665) and the ready to use phospho specific monoclonal antibody labeled with Europium Cryptate (Eu(K)) are kept constant. The optimal enzyme concentration is chosen at EC_{80} of the titration curve obtained

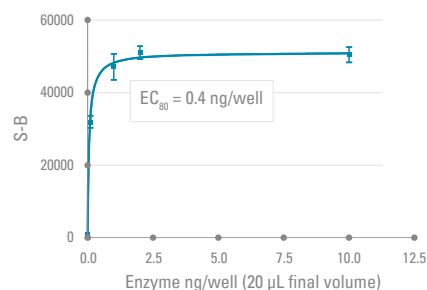


Fig. 2: MAPKAP-K2 titration

2- ENZYME KINETIC

Enzyme kinetic depends on the kinase and the substrate concentrations. A time course study is performed using 0.4 ng of MAPKAP-K2 per well, the constant concentration of kinase determined in the previous experiment, 1 μ M of substrate and a non-limiting ATP concentration (100 μ M). The reaction is stopped at different end points by the addition of the detection reagents (at 1, 2, 5, 10, 15, 30 and 60 min). The optimal incubation period for MAPKAP-K2 (0.4 ng/well) to achieve maximum signal and a linear time course is chosen at 5 min, this incubation time is kept constant for the rest of the optimization.

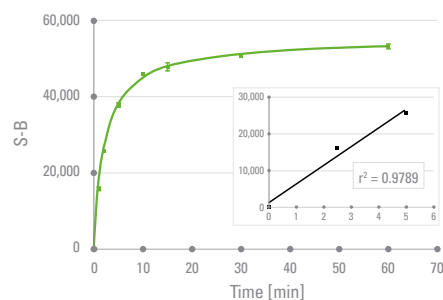


Fig 3: MAPKAP-K2 enzyme kinetic

3- SUBSTRATE TITRATION

In order to determine the substrate K_M (app), assays are run under the conditions previously obtained (enzyme concentration: 0.4 ng/well and incubation period: 5 min) using substrate concentrations ranging from 1 nM to 2 μ M. During the detection step, the SA-XL665 concentration is adjusted to keep the substrate/streptavidin-XL665 ratio constant at 8/1. The signal is plotted versus substrate concentrations, and the K_M (app) of 80 nM is calculated using Michaelis-Menten equation.

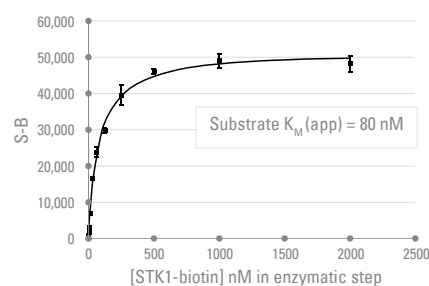


Fig 4: MAPKAP-K2 substrate titration.

4- ATP TITRATION

Assays are run at a non-limiting substrate concentration with ATP from 1.7 nM to 300 μ M, while the enzyme quantity (0.4 ng/well) and the incubation period (5 min) are kept constant.

As in the previous step, the ATP K_M (app) of 2.3 μ M is calculated from the resulting plot of the signal versus ATP concentrations.

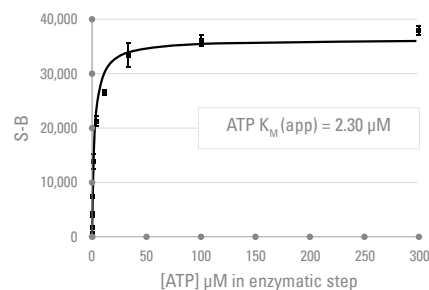


Fig 5: MAPKAP-K2 ATP titration

5- DETECTION STEP OPTIMIZATION

The optimization of substrate-biotin/SA-XL665 ratio is an important step which may lead to a substantial increase in signal. The assay is performed using the optimal enzyme, ATP and substrate concentrations. Three different molar ratios of substrate-biotin/streptavidin-XL665 are tested (2/1, 4/1, 8/1). The optimal ratio obtains a good compromise between signal level and reagent consumption.

6- INHIBITOR TITRATION

The kinase activity is tested over a broad range of inhibitor concentrations to generate a dose-response curve. The test is generally run using the optimal assay conditions as determined previously. Table beside shows reference inhibitor IC_{50} for STK and TK kinases using HTRF KinEASE.

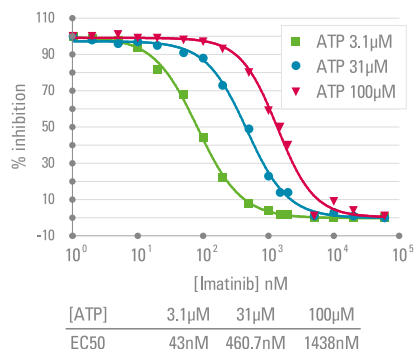


Fig 6: Imatinib is an ATP dependant PKI. The imatinib IC_{50} is calculated with HTRF KinEASE at increasing ATP concentrations, K_M (3.1 μM), 10 K_M (31 μM), an excess of ATP (100 μM), 1 μM TK substrate-biotin, 1 ng/well enzyme, 4/1 ratio biotin/SaXL665, [SEB]=5 nM(1) and 20 min incubation for the enzymatic step. The results obtained are consistent with pharmacological data already published.

ENZYME	HTRF KINEASE	COMPOUND	IC_{50} NM
Abl	TK	Staurosporine Imatinib*	108 43*
Csk	TK	Staurosporine Sorafenib*	493 4,220
EGFR	TK	Staurosporine Gefitinib*	1,07 1*
JAK2	TK	Staurosporine Sunitinib*	0.30 800
JAK3	TK	Staurosporine Sunitinib*	1 7,64
PTK5	TK	Staurosporine Sorafenib*	39 >100
MAPKAP-K2	STK S1	Staurosporine	92
CaMK IV	STK S1	Staurosporine	4.3
PKC beta 2	STK S1	Staurosporine	6
Rsk3	STK S1	Staurosporine	3.1
Rock-II	STK S2	Staurosporine	10.4
Pim-1	STK S2	Staurosporine	8.7

Table 1: The IC_{50} of TKIs of interest are calculated with HTRF KinEASE on a selection of kinases. The values obtained are similar to the ^{32}P incorporation method used as a reference.

(1) SEB (Cisbio proprietary buffer) can be used for optimal kinase TK activity and is provided with KinEASE TK.

A PLATFORM VALIDATED ON MORE THAN 160 KINASES

Today, the platform has been validated on more than 100 Ser/Thr kinases and 60 Tyrosine kinases including both receptor and cytoplasmic kinases. Please refer to www.htrf.com/kaa for an updated list of kinases already tested with HTRF. Serine threonine kinases validated with HTRF KinEASE STK & TK:

HTRF® KinEASE STK S1	AMPK	BrSK2	CAMK1	CaMKII α	CaMKII β	CaMKII δ	CaMKII γ	CaMKIV	CHEK1	CHEK2	CHK1
	CHK2	DAPK1	DAPK2	DAPK3	DCAMKL2	DRAK1	eEF-2K	LKB1	MAPKAP-K1 α	MAPKAPK1 β	MAPKAP-K2
	MAPKAP-K3	MAPKAPK5	MELK	MLCK	Mnk2	PASK	PHK γ 2	PKC α	PKC β I	PKC β II	PKC δ
	PKC ε	PKC γ	PKC ι	PKC μ	PKC θ	PKC ζ	PKD2	PRAK	PRK2	PRKA	PRKD2
	PRKG2	RPS6KA2	Rsk1	Rsk2	Rsk3	STK22B	STK22D	TSSK1	TSSK2	ZIPK	
HTRF® KinEASE STK S2	Aurora A	Aurora B	MARK2	NEK11	PAK2	PAK3	PAK4	PAK5	PAK6	PAR-1B α	PKA
	PKG1 α	PKG1 β	PRKACA	ROCK-I	ROCK-II	ROK α	RPS6KA6	Rsk4	STK6		
HTRF® KinEASE STK S3	AKT1	AKT2	AKT3	ARK5	ASK1	Aurora C	BrSK1	CDC42 BPA	CDC42 BPB	CLK3	COT
	DMPK	DYRK2	DRAK2	GRK5	GRK6	HIPK2	HIPK3	IKK α	IKK β	LOK	MAP3K8
	MARK1	MINK	MINK1	MRCK α	MRCK β	MSK1	MSK2	MSSK1	MST1	MST2	NEK2
	NEK3	NEK6	NEK7	NLK	p70S6K	Pim-1	Pim-2	PKB α	PKB β	PKB γ	PLK3
	RPS6KA4	RPS6KA5	RPS6KB1	SAD1	SGK1	SGK2	SGK3	SGKL	SIK	Snk	STK23
	STK3	STK4	TBK1	WNK2	WNK3						
HTRF® KinEASE TK	Abl	ABL2	ALK	Arg	Axl	Blk	Bmx	BRK	BTK	c-Kit	CSF1R
	CSK	c-SRC	DDR2	EGFR	EphA2	EphA3	EphA4	EphA5	EphA7	EphA8	EphB1
	EphB2	EphB3	EphB4	ErbB4	FAK	Fer	Fes	FGFR1	FGFR2	FGFR3	FGFR4
	Fgr	Flt1	Flt3	Flt4	Fms	Fps	FRK	Fyn	Hck	IGF-1R	INSRR
	Insulin R	IRR	ITK	JAK2	JAK3	KDR	Kit	LCK	Lyn	Mer	Met
	MST1R	MuSK	NTRK1	NTRK2	PDGFR α	PDGFR β	PTK2	PYK2	PTK5	PTK6	Ret
	Ron	Ros	Rse	Src	Syk	TEK	Tie2	TRKA	TRKB	TYRO3	VEGFR1
	VEGFR2	VEGFR3	Yes	YES1	ZAP-70						

The recommended substrates were determined for Millipore kinases.

CONCLUSION

Cisbio's line of HTRF KinEASE kits is based on our patented HTRF technology and can be used as a universal tool for assessing Serine/Threonine and Tyrosine kinase activity.

HTRF KinEASE kits limit assay development time and are easily miniaturizable and flexible, meaning the assay can be performed under a wide range of kinase assay conditions, for instance with low consumption of enzyme or with any ATP concentration.

RELATED INFORMATION

HTRF® KinEASE™ TK: A new solution for Tyrosine kinases screening.

Drexler C., SBS 13th annual conference 2007, Montreal (Canada).

Universal Expanded Platform to address serine/threonine and Tyrosine Kinases.

Tardieu J.L., BIForum Europe, 42 (6/2007).

HTRF® KinEASE™ : a new solution for screening serine-threonine kinases.

Drexler C., Nature Methods, June 2006.

HTRF® KinEASE™: Development of sensitive, reliable Aurora A and AMPK kinase inhibitory assays.

Martinez S., SBS 12th annual conference. September 2006, Seattle (USA).

HTRF® KinEASE™: A universal assay for Serine/Threonine kinases.

Claret E., SBS 12th annual conference. September 2006, Seattle (USA).

ORDERING INFORMATION

HTRF® KinEASE™ for Serine / Threonine and Tyrosine kinases

DESCRIPTION	TESTS	PART#
HTRF® KinEASE™-STK discovery (STK substrates 1, 2 and 3-biotin)	1,000 tests	62ST0PEB
HTRF® KinEASE™-STK S1 (STK substrate 1-biotin)	1,000 tests	62ST1PEB
	20,000 tests	62ST1PEC
	100,000 tests	62ST1PEJ
HTRF® KinEASE™-STK S2 (STK substrate 2-biotin)	1,000 tests	62ST2PEB
	20,000 tests	62ST2PEC
	100,000 tests	62ST2PEJ
HTRF® KinEASE™-STK S3 (STK substrate 3-biotin)	1,000 tests	62ST3PEB
	20,000 tests	62ST3PEC
	100,000 tests	62ST3PEJ
HTRF® KinEASE™-TK (TK substrate biotin)	1,000 tests	62TK0PEB
	20,000 tests	62TK0PEC
	100,000 tests	62TK0PEJ

HTRF® KinEASE™ companion component

DESCRIPTION	TESTS	PART#
STK substrate 1-biotin	50 µg/vial	61ST1BLE
	500 µg/vial	61ST1BLC
STK substrate 2-biotin	50 µg/vial	61ST2BLE
	500 µg/vial	61ST2BLC
STK substrate 3-biotin	50 µg/vial	61ST3BLE
	500 µg/vial	61ST3BLC
TK substrate -biotin	50 µg/vial	61TK0BLE
	500 µg/vial	61TK0BLC
Sa-XL665	250 µg	610SAXLA
	1 mg	610SAXLB
	3 mg	610SAXLG
5x Enzymatic buffer	50 mL	62EZBFDD
SEB buffer for HTRF KinEASE TK assays	20,000 tests (lyoph, to be reconst. with 5 ml H2O)	61SEBALB
HTRF® Detection buffer	200 mL	62SDBRDF

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