

HTRF® Research Reagents

HTRF KinEASE STK S1 kit - 1K pts

Product Number : 62ST1PEB Lot Number : 11K

Expiration date : 2026-04-18 Document version : 1

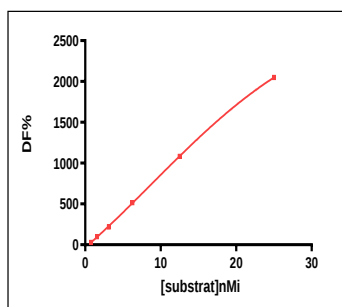
Reagents :	Lot n°
HTRF STK-Substrate 1-biotin - 50 µg	08RA
HTRF Streptavidin-XL665 - 5K pts	183A
HTRF Enzymatic Buf./kinase (5X) - 10 mL	16AR
HTRF KinEASE STK kit - 1K pts Eu	11RA
HTRF KinEASE Detect. Buf. - 40 mL	20A
/	/
/	/

Results :

Incubation : 1h at RT
Plate : Costar 96 wells half volume black
Reader : Pherastar Plus

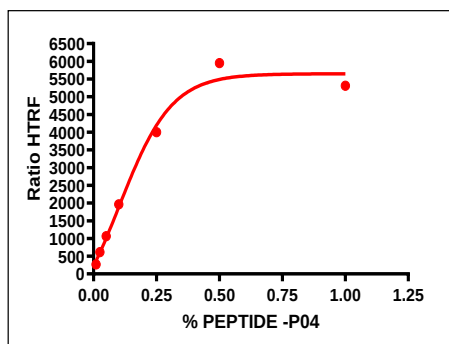
Results : Substrate validation

[S1-biotin]nMi	DF%	Specifications Rev 02
0,781	29%	
1,563	98%	
3,125	218%	
6,25	515%	
12,5	1083%	
25	2048%	≥ 1200
Ratio negative 25nM	369	≤ 600
Ratio cryptate blank	347	≤ 600
EC 50	12,2	11 – 14 nMi



Results : Cryptate and XL validation

% Peptide PO4	DF (%)	Specifications Rev 02
0,01	269%	
0,025	613%	
0,05	1065%	
0,1	1963%	
0,25	3999%	
0,5	5948%	
1	5309%	≥ 3500
Ratio negative	468	≤ 700
Ratio Cryptate Blank	438	≤ 600
EC 50	0,25	0.15 - 0.33 nMi



HTRF KinEASE STK S1 kit - 1K pts - Lot n° : 11K

☒ Accepted☐ Rejected

Quality control date : 2025-06-16 Supervisor's initials : MP

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.