

HTRF® Research Reagents

HTRF KinEASE STK S2 kit - 1K pts

Product Number : 62ST2PEB Lot Number : 11H

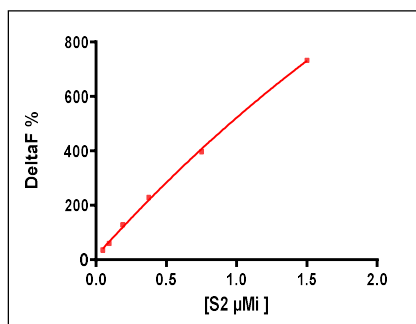
Expiration date : 2026-04-18 Document version : 1

Reagents :	Lot n°
HTRF STK-Substrate 2-biotin - 50 µg	11A
HTRF Streptavidin-XL665 - 5K pts	181AR
HTRF Enzymatic Buf./kinase (5X) - 10 mL	16AR
HTRF KinEASE STK kit - 1K pts Eu	11RA
HTRF KinEASE Detect. Buf. - 40 mL	18RA

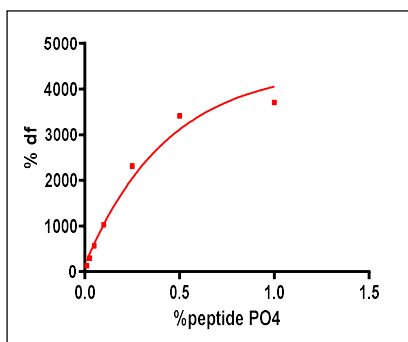
Results :

Incubation : 1h at RT
Plate : Costar 96 wells half volume black
Reader : Envision
Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

[S2-biotin]µMi	DF%	Specifications Rev 02
0,047	179%	
0,094	343%	
0,188	612%	
0,375	1001%	
0,75	1664%	
1,5	1808%	≥ 700
Ratio negative 1.5µM	765	≤ 800
Ratio Cryptate Blank	403	≤ 600
EC50	0,45	0.35 – 0.7 µMi



% Peptide PO4	DF (%)	Specifications Rev 02
0,01	214%	
0,025	460%	
0,05	774%	
0,1	1290%	
0,25	2746%	
0,5	3793%	
1	4087%	≥3500
	20,4354579	
Ration negative	432	≤700
Ratio Cryptate Blank	382	≤ 600
EC 50	0,17	0.15 -0.33 nMi



HTRF KinEASE STK S2 kit - 1K pts - Lot n° : 11H

☒ Accepted☐ Rejected

Quality control date : 2024-07-02 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.