

HTRF® Research Reagents

HTRF KinEASE STK S3 kit - 1K pts

Product Number : 62ST3PEB Lot Number : 11EA

Expiration date : 2025-02-02 Document version : 1

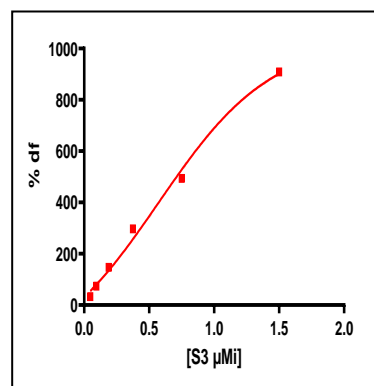
Reagents :	Lot n°
HTRF STK-Substrate 3-biotin - 50 µg	08RA
HTRF Streptavidin-XL665 - 5K pts	181AR
HTRF Enzymatic Buf./kinase (5X) - 10 mL	14RA
HTRF KinEASE STK kit - 1K pts Eu	11RA
HTRF KinEASE Detect. Buf. - 40 mL	17RA

Results :

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

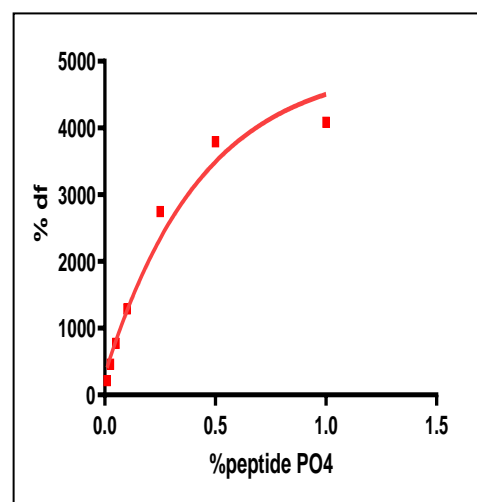
Results : Substrate validation

[S3-biotin]µMi	DF%	Specifications Rev 02
0,047	33%	
0,094	74%	
0,188	147%	
0,375	297%	
0,75	494%	
1,5	909%	≥ 700
Ratio negative 1.5µM	498	≤1000
Ratio cryptate blank	370	≤600
EC 50	0,65	0.35 – 0.7 µMi



Results : Cryptate and XL validation

% 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
Ration negative	432	≤700
Ratio Cryptate Blank	382	≤ 600
EC 50	0,17	0.15 -0.33 nMi



HTRF KinEASE STK S3 kit - 1K pts - Lot n° : 11EA

Accepted

Rejected

Quality control date : 2023-12-08 Supervisor's initials : PR

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.