

HTRF® Research Reagents

HTRF KinEASE STK S3 kit - 20K pts

Product Number : 62ST3PEC Lot Number : 25CA

Expiration date : 2025-08-03 Document version : 1

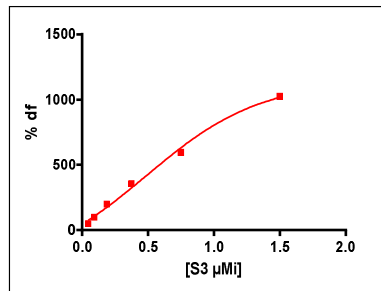
Reagents :	Lot n°
HTRF STK-Substrate 3-biotin - 500 µg	14RA
HTRF Streptavidin-XL665 - 3 mg	181AR
HTRF KinEASE STK kit - 20K pts Eu	25RA
HTRF Enzymatic Buf. kinase (5X) - 50 mL	15RA
HTRF KinEASE Detect. Buf. - 200 mL	19RA

Results :

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

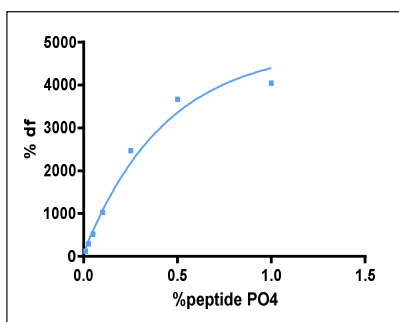
Results : Substrate validation

[S3-biotine] µMi	DF%	Specification Rev 02
0.047	49%	
0.094	98%	
0.188	198%	
0.375	355%	
0.75	595%	
1.5	1026%	≥ 700
Ratio negative 1.5µM	483	≤ 1000
Ratio cryptate blank	362	≤ 600
EC 50	0.6	0.35 - 0.7 µMi



Results : Cryptate and XL validation

% Peptide PO4	DF (%)	Specification Rev 02
0.01	125%	
0.025	294%	
0.05	524%	
0.1	1032%	
0.25	2464%	
0.5	3668%	
1	4046%	≥ 3500
Ratio negative	444	≤ 700
Ratio Cryptate Blank	450	≤ 600
EC 50	0.23	0.15 - 0.33 nMi



HTRF KinEASE STK S3 kit - 20K pts - Lot n° : 25CA

Accepted

Rejected

Quality control date : 2023-10-28	Supervisor's initials : FL
-----------------------------------	----------------------------

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.