

HTRF® Research Reagents

HTRF KinEASE STK Discovery kit - 1K pts

Product Number : 62ST0PEB Lot Number : 11H

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

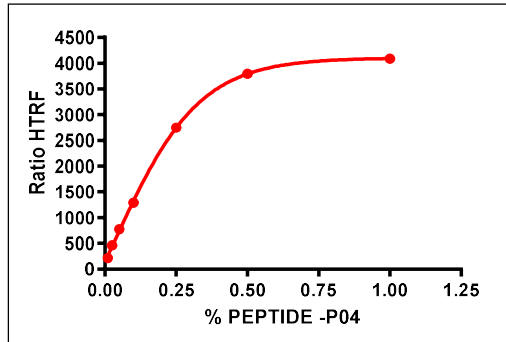
Reagents :	Lot n°
HTRF STK-Substrate 1-biotin - 50 µg	08RA
HTRF STK-Substrate 2-biotin - 50 µg	11A
HTRF STK-Substrate 3-biotin - 50 µg	08RA
HTRF Streptavidin-XL665 - 5K pts	181AR
HTRF Enzymatic Buf./kinase (5X) - 10 mL	16AR
HTRF KinEASE STK kit - 1K pts Eu	11RA
HTRF KinEASE Detect. Buf. - 40 mL	17RA

Results :

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

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



Assay characteristics :

	Found value	Specifications Rev 01
DF % max	4087	≥3500
Ratio negative	432	≤700
Ratio cryptate Blanc	382	≤ 600
EC50	0,17	0.15 -0.33 nMi

HTRF KinEASE STK Discovery 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.