

HTRF® Research Reagents

HTRF KinEASE TK kit - 1K pts

Product Number : 62TK0PEB Lot Number : 09E

Expiration date : 2025-10-01 Document version : 1

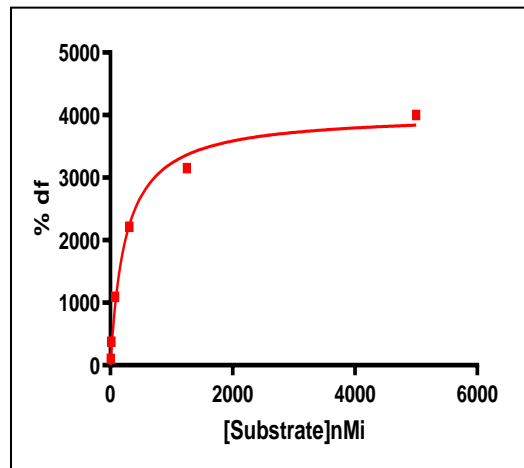
Reagents :	Lot n°
HTRF Streptavidin-XL665 - 5K pts	182A
HTRF Enzymatic Buf./kinase (5X) - 10 mL	16AR
HTRF KinEASE Detect. Buf. - 40 mL	17RA
HTRF Suppl. Enzymatic Buf. - 2.5 nmoles	12RA
HTRF KinEASE TK kit - 1K pts Eu	09AR
HTRF TK Substrate-biotin - 50 µg	09RA

Results :

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

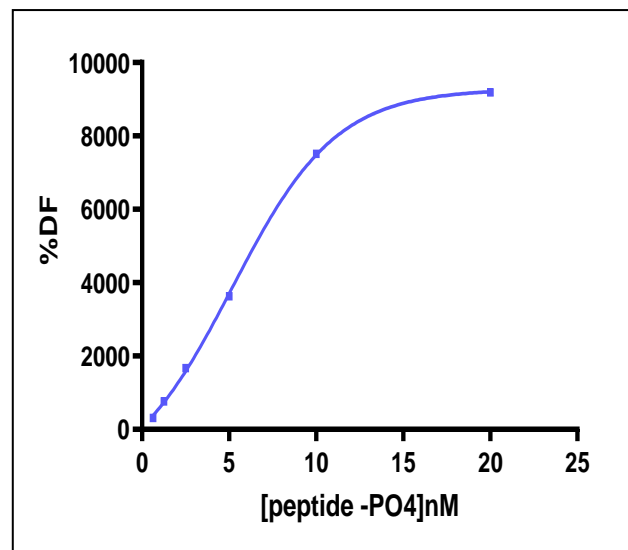
Results : Substrate validation

[TK Sub-biotin] nMi	DF (%)	Specifications Rev 03
4,88	105%	
19,53	378%	
78,1	1095%	
312,5	2215%	
1250	3153%	
5000	4003%	≥ 3500
Ratio negative 5000 nMi	515	≤ 1000
Ratio Cryptate Blank	339	≤ 600



Results : Cypate and XL validation

% peptide PO4	DF %	Specifications Rev 03
0,625	312%	
1,25	759%	
2,5	1675%	
5	3627%	
10	7511%	
20	9190%	≥ 5000%
Ratio negative	376	≤ 600
Ratio Cryptate Blank	337	≤ 600
EC 50	5,9	4 – 9 nMi



HTRF KinEASE TK kit - 1K pts - Lot n° : 09E

Accepted Rejected

Quality control date : 2025-01-10	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.