

HTRF® Research Reagents

HTRF KinEASE TK kit - 1K pts

Product Number : 62TK0PEB Lot Number : 08K

Expiration date : 2024-03-31 Document version : 1

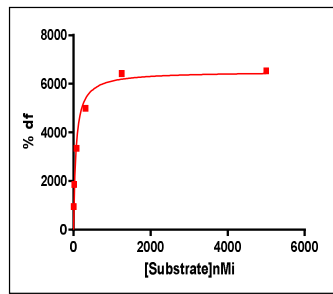
Reagents :	Lot n°
HTRF Streptavidin-XL665 - 5K pts	175A
HTRF Enzymatic Buf./kinase (5X) - 10 mL	14RA
HTRF KinEASE Detect. Buf. - 40 mL	17 A
HTRF Suppl. Enzymatic Buf. - 2.5 nmoles	12A
HTRF KinEASE TK kit - 1K pts Eu	08A
HTRF TK Substrate-biotin - 50 µg	08A
HTRF KinEASE TK kit - modop	5

Results :

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

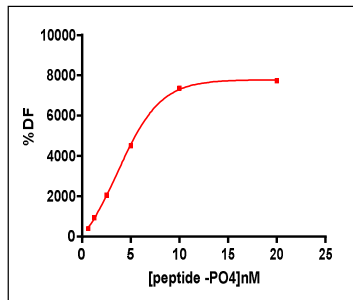
Results : Substrate validation

[TK Sub-biotin] nMi	DF (%)	Specifications Rev 03
4.88	958%	
19.53	1857%	
78.1	3348%	
312.5	4987%	
1250	6421%	
5000	6536%	≥ 3500
Ratio negative 5000 nMi	605	≤ 1000
Ratio Cryptate Blank	351	≤ 600



Results : Cypate and XL validation

% peptide PO4	DF %	Specifications Rev 03
0.625	401%	
1.25	940%	
2.5	2063%	
5	4516%	
10	7367%	
20	7739%	≥ 5000%
Ratio negative	404	≤ 600
Ratio Cryptate Blank	380	≤ 600
EC 50	4.3	4 – 9 nMi



HTRF KinEASE TK kit - 1K pts - Lot n° : 08K

Accepted  Rejected

Quality control date : 2023-09-06	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.