

HTRF® Research Reagents

HTRF KinEASE TK kit - 20K pts

Product Number : 62TK0PEC Lot Number : 16B

Expiration date : 2027-01-08 Document version : 1

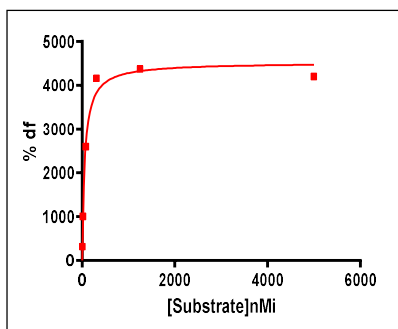
Reagents :		Lot n°
HTRF Streptavidin-XL665 - 3 mg		182A
HTRF Enzymatic Buf. kinase (5X) - 50 mL		15RA
HTRF KinEASE Detect. Buf. - 200 mL		19RA
HTRF Suppl. Enzymatic Buf. - 12.5 nmoles		11A
HTRF KinEASE TK kit - 20K pts Eu		16A
HTRF TK Substrate-biotin - 500 µg		24A

Results :

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

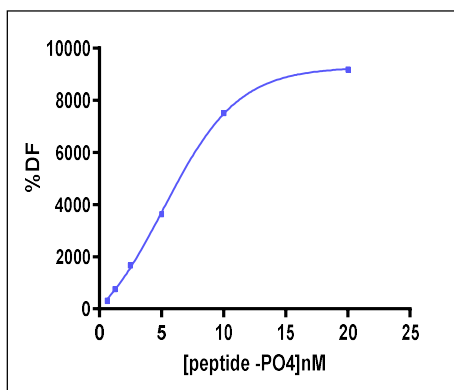
Results : Substrate validation

[TK Sub-biotin] nMi	DF (%)	Specifications Rev 03
4,88	313%	
19,53	1007%	
78,1	2603%	
312,5	4165%	
1250	4379%	
5000	4202%	≥ 3500
Ratio negative 5000 nMi	709	≤ 1000
Ratio Cryptate Blank	414	≤ 600



Results : Cypate and XL validation

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



HTRF KinEASE TK kit - 20K pts - Lot n° : 16B

Accepted Rejected

Quality control date : 2024-09-09	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.