

Quality Control report

HTRF KinEASE TK kit - 1K pts
 (62TKOPEB)
 Lot n°08J
 Expiration date : 2024-03-31

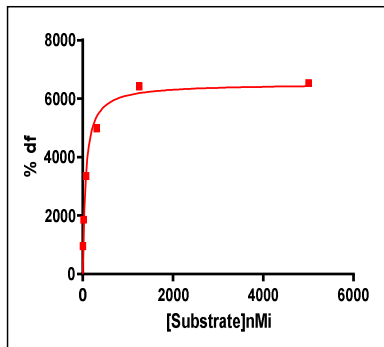
Reagents :	Lot n°
HTRF Streptavidin-XL665 - 5K pts	175A
HTRF Enzymatic Buf./kinase (5X) - 10 mL	14A
HTRF KinEASE Detect. Buf. - 40 mL	17 A
HTRF Suppl. Enzymatic Buf. - 2.5 nmoles	12A
HTRF KinEASE TK kit - 1K pts Eu	08A
HTRF TK Substrate-biotin - 50 µg	08A
HTRF KinEASE TK kit - modop	5

Results :

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

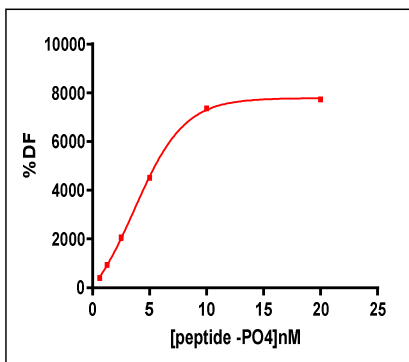
Results : Substrate validation

[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°08J

Accepted

Rejected

Quality control date : 2023-07-13 Supervisor's initials : FL

