

## HTRF® Research Reagents

## HTRF KinEASE TK kit - 1K pts

Product Number : 62TK0PEB Lot Number : 09H

Expiration date : 2026-11-08 Document version : 1

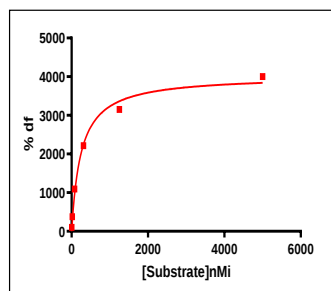
| Reagents :                              | Lot n° |
|-----------------------------------------|--------|
| HTRF Streptavidin-XL665 - 5K pts        | 183A   |
| HTRF Enzymatic Buf./kinase (5X) - 10 mL | 16AR   |
| HTRF KinEASE Detect. Buf. - 40 mL       | 20A    |
| HTRF Suppl. Enzymatic Buf. - 2.5 nmoles | 12RA   |
| HTRF KinEASE TK kit - 1K pts Eu         | 09AR   |
| HTRF TK Substrate-biotin - 50 µg        | 09RA   |
| /                                       | /      |
| /                                       | /      |

## Results :

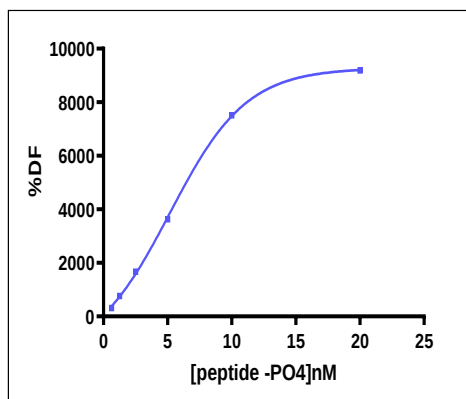
Incubation : 1h at RT  
 Plate : 96 wells half volume black  
 Reader : Pherastar Plus

## Results : Substrate validation

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

Results : Cryptate  
and XL  
validation

| % peptide PO4           | DF %  | Specifications<br>Rev 03 |
|-------------------------|-------|--------------------------|
| 0,625                   | 312%  |                          |
| 1,25                    | 759%  |                          |
| 2,5                     | 1675% |                          |
| 5                       | 3627% |                          |
| 10                      | 7511% |                          |
| 20                      | 9190% | ≥ 5000%                  |
| Ratio negative          | 539   | ≤ 600                    |
| Ratio Cryptate<br>Blank | 455   | ≤ 600                    |
| EC 50                   | 5,6   | 4 – 9 nMi                |



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

☒ Accepted☐ Rejected

Quality control date : 2025-07-08 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.