

HTRF® Research Reagents

HTRF PD1/PD-L1 Bind. kit - 500 pts

Product Number : 64PD1PEG Lot Number : 07E

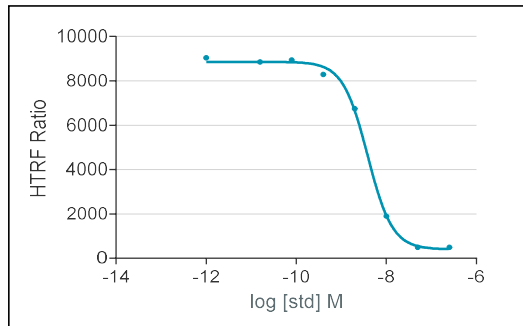
Expiration date : 2026-04-24 Document version : 1

| Reagents : | Lot n° |
|--|--------|
| HTRF PPI-Europium Detect. Buf. - 20 mL | 08A |
| HTRF PD1/PD-L1 Bind. kit - 500 pts Tag2-PD1 | 06RA |
| HTRF PD1/PD-L1 Bind. kit - Std | 05A |
| HTRF PD1/PD-L1 Bind. kit - 500 pts Eu Cryptate reagent | 07A |
| HTRF PD1/PD-L1 Bind. kit - 500 pts Tag1-PD-L1 | 06RA |
| HTRF PD1/PD-L1 Bind. kit - 500 pts XL665 antibody | 07A |

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

| PD1/PD- L1 | Conc. (nMf) | RATIO |
|------------|-------------|-------|
| Std 0 | 0 | 9044 |
| Std 1 | 0,016 | 8854 |
| Std 2 | 0,08 | 8943 |
| Std 3 | 0,4 | 8299 |
| Std 4 | 2 | 6747 |
| Std 5 | 10 | 1901 |
| Std 6 | 50 | 494 |
| Std 7 | 250 | 499 |



Assay characteristics :

| | Found value | Specifications Rev 01 |
|-------|-------------|-----------------------|
| IC 50 | 2,87 | 2 - 7 nM |
| S/B | 18 | 16-26 |

HTRF PD1/PD-L1 Bind. kit - 500 pts - Lot n° : 07E

Accepted

Rejected

| | |
|-----------------------------------|----------------------------|
| Quality control date : 2024-11-12 | 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.