

HTRF® Research Reagents

HTRF IL23/IL23R Bind. kit - 500 pts

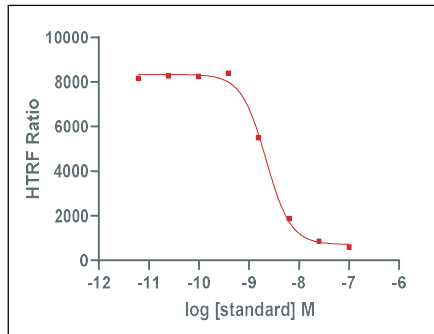
Product Number : 64BDPIL23PEG Lot Number : 04B

Expiration date : 2026-02-07 Document version : 1

| Reagents : | Lot n° |
|--|--------|
| HTRF PPI-Europium Detect. Buf. - 20 mL | 07RA |
| HTRF IL23/IL23R Bind. kit - 500 pts Tag2-IL23R | 05RA |
| HTRF IL23/IL23R Bind. kit - Std | 03A |
| HTRF IL23/IL23R Bind. kit - 500 pts d2 reagent | 04A |
| HTRF IL23/IL23R Bind. kit - 500 pts Eu Cryptate antibody | 04A |
| HTRF IL23/IL23R Bind. kit - 500 pts Tag1-IL23 | 05RA |
| | |
| | |

Results : 2h at RT
 Incubation : 384 wells small volume white plate 20 µL
 Plate : Pherastar Plus
 Reader : log(inhibitor) vs. response -- Variable slope (four parameters)
 Fitting model :

| std | Conc. (pM) | RATIO |
|-------|------------|-------|
| Std 0 | 0 | 8 161 |
| Std 1 | 24,4 | 8 268 |
| Std 2 | 97,7 | 8 247 |
| Std 3 | 390 | 8 384 |
| Std 4 | 1563 | 5 512 |
| Std 5 | 6250 | 1 877 |
| Std 6 | 25000 | 861 |
| Std 7 | 100000 | 598 |



Assay characteristics :

| | Found value | Specifications Rev 01 |
|---------|-------------|-----------------------|
| negatif | 416 | / |
| S/B | 20 | 15-30 |
| IC50 | 2,19 | 1.5 - 3.4 nM |

HTRF IL23/IL23R Bind. kit - 500 pts - Lot n° : 04B

Accepted

Rejected

| | | | |
|------------------------|------------|-------------------------|----|
| Quality control date : | 2024-06-25 | 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.