

HTRF® Research Reagents

HTRF Cereblon Bind. kit - 10K pts

Product Number : 64BDCRBNPEH Lot Number : 06A

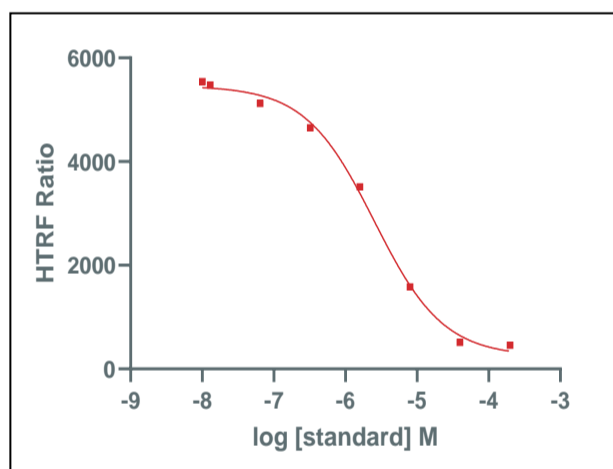
Expiration date : 2025-10-03 Document version : 1

| Reagents :   | Lot n° |
|--|--------|
| HTRF Diluent 9 (5X) - 100 mL   | 04RA   |
| HTRF PROTAC Bind. Buf. 1 - 200 mL                                      | 04RA   |
| HTRF Cereblon Bind. kit - 10K pts Eu Cryptate antibody                 | 06RA   |
| HTRF Cereblon Bind. kit - 10K pts Human cereblon WT GST-tagged protein | 11A    |
| HTRF Cereblon Bind. kit - 10K pts Thalidomide Red                      | 06A    |
| HTRF Cereblon Bind. kit - Std  | 04RA   |

Results :

Incubation : 3h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : log(inhibitor) vs. response -- Variable slope (four parameters)

| std   | Conc. (µM) | RATIO |
|-------|------------|-------|
| Std 0 | 0          | 5543  |
| Std 1 | 0,0128     | 5478  |
| Std 2 | 0,064      | 5123  |
| Std 3 | 0,32       | 4652  |
| Std 4 | 1,6        | 3511  |
| Std 5 | 8          | 1583  |
| Std 6 | 40         | 511   |
| Std 7 | 200        | 457   |



Assay characteristics :

|      | Found value | Specifications Rev 01 |
|------|-------------|-----------------------|
| S/B  | 12          | 12 -18                |
| IC50 | 2,5         | 1 - 4 µM              |

HTRF Cereblon Bind. kit - 10K pts - Lot n° : 06A

Accepted

Rejected

|                                   |                            |
|-----------------------------------|----------------------------|
| Quality control date : 2024-02-23 | Supervisor's initials : FL |
|-----------------------------------|----------------------------|

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.