

HTRF® Research Reagents

HTRF Cereblon Bind. kit - 10K pts

Product Number : 64BDCRBNEPH Lot Number : 04F

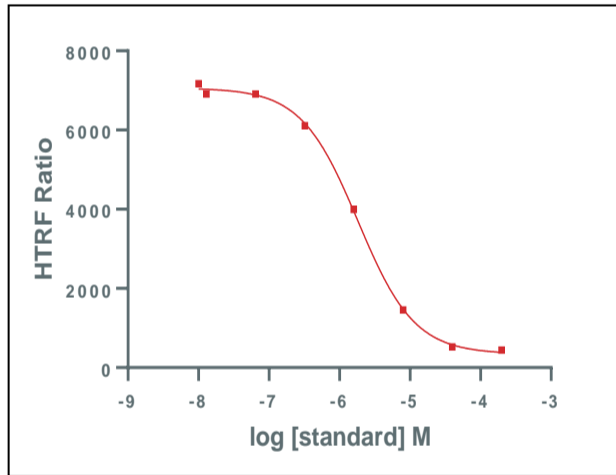
Expiration date : 2025-01-08 Document version : 1

Reagents :	Lot n°
HTRF Diluent 9 (5X) - 100 mL	03RA
HTRF PROTAC Bind. Buf. 1 - 200 mL	04RA
HTRF Cereblon Bind. kit - 10K pts Eu Cryptate antibody	03RA
HTRF Cereblon Bind. kit - 10K pts Human cereblon WT GST-tagged protein	11A
HTRF Cereblon Bind. kit - 10K pts Thalidomide Red	04RA
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	5548
Std 1	0,0128	5219
Std 2	0,064	5137
Std 3	0,32	4685
Std 4	1,6	3284
Std 5	8	1293
Std 6	40	513
Std 7	200	467



Assay characteristics :

	Found value	Specifications Rev 01
S/B	12	12 -18
IC50	2,15	1 - 4 µM

HTRF Cereblon Bind. kit - 10K pts - Lot n° : 04F

Accepted

Rejected

Quality control date : 2024-02-08	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.