

HTRF® Research Reagents

HTRF Cereblon Bind. kit - 10K pts

Product Number : 64BDCRBNEPH Lot Number : 06D

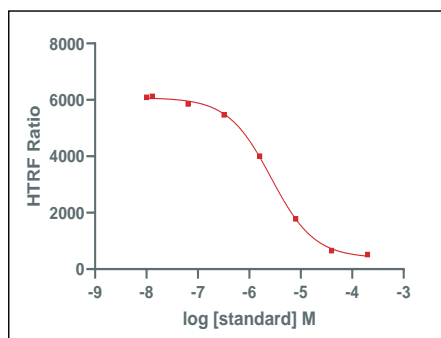
Expiration date : 2026-03-28 Document version : 1

Reagents :	Lot n°
HTRF Diluent 9 (5X) - 100 mL	04RA
HTRF PROTAC Bind. Buf. 1 - 200 mL	05AR
HTRF Cereblon Bind. kit - 10K pts Eu Cryptate antibody	06RA
HTRF Cereblon Bind. kit - 10K pts Human cereblon WT GST-tagged protein	14A
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	6089
Std 1	0,0128	6124
Std 2	0,064	5856
Std 3	0,32	5473
Std 4	1,6	4002
Std 5	8	1784
Std 6	40	658
Std 7	200	518



Assay characteristics :

	Found value	Specifications Rev 01
S/B	12	12 -18
IC50	2,6	1 - 4 µM

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

Accepted

Rejected

Quality control date :	2025-03-21	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.