

HTRF® Research Reagents

HTRF Cereblon Bind. kit - 500 pts

Product Number : 64BDCRBNPEG Lot Number : 07F

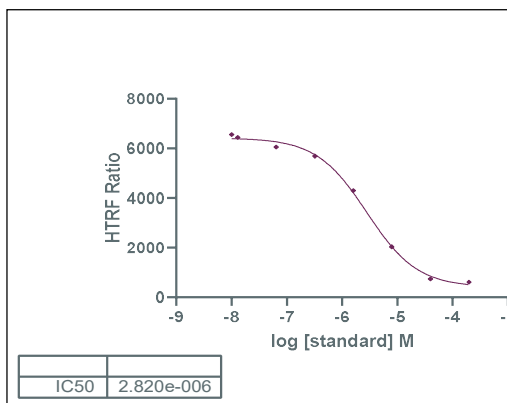
Expiration date : 2025-10-03 Document version : 1

Reagents :		Lot n°
HTRF Diluent 9 (5X) - 2 mL		04A
HTRF PROTAC Binding Buffer 1 - 20 mL		05AR
HTRF Cereblon Bind. kit - Std		04RA
HTRF Cereblon Bind. kit - 500 pts Eu Cryptate antibody		07AR
HTRF Cereblon Bind. kit - 500 pts Human cereblon WT GST-tagged protein		12A
HTRF Cereblon Bind. kit - 500 pts Thalidomide Red		07BR

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	6559
Std 1	0,0128	6441
Std 2	0,064	6056
Std 3	0,32	5688
Std 4	1,6	4302
Std 5	8	2034
Std 6	40	740
Std 7	200	612



Assay characteristics :

	Found value	Specifications Rev 01
S/B	11	
IC50	2,8	1 - 4 µM

HTRF Cereblon Bind. kit - 500 pts - Lot n° : 07F

☒ Accepted☐ Rejected

Quality control date :	2024-10-05	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.