

HTRF® Research Reagents

HTRF Cereblon Bind. kit - 500 pts

Product Number : 64BDCRBNPEG Lot Number : 08B

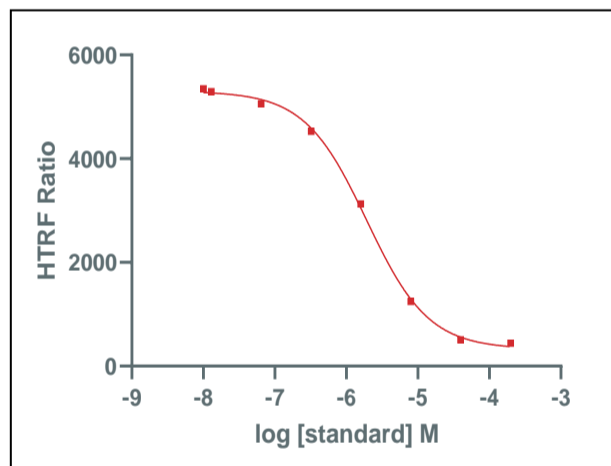
Expiration date : 2025-01-08 Document version : 1

Reagents :	Lot n°
HTRF Diluent 9 (5X) - 2 mL	03B
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	09RA
HTRF Cereblon Bind. kit - 500 pts Thalidomide Red	08A

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	5455
Std 1	0,0128	5423
Std 2	0,064	5161
Std 3	0,32	4799
Std 4	1,6	3457
Std 5	8	1671
Std 6	40	534
Std 7	200	450



Assay characteristics :

	Found value	Specifications Rev 01
S/B	12	12 -18
IC50	2,64	1 - 4 µM

HTRF Cereblon Bind. kit - 500 pts - Lot n° : 08B

Accepted

Rejected

Quality control date : 2024-01-03	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.