

HTRF® Research Reagents

HTRF MDM2 Binding Kit - 500 pts

Product Number : 64BDMDM2PEG Lot Number : 01E

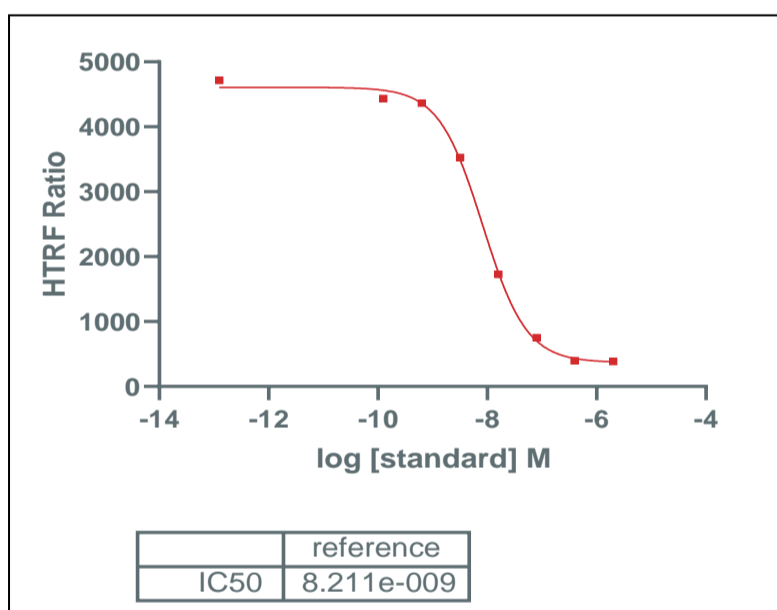
Expiration date : 2026-07-01 Document version : 1

Reagents :	Lot n°
HTRF Diluent 9 (5X) - 2 mL	04A
HTRF PROTAC Binding Buffer 3 - 20 mL	01RA
HTRF MDM2 Binding Kit - Std	01RA
HTRF MDM2 Binding Kit - 500 pts Eu Cryptate antibody	01RA
HTRF MDM2 Binding Kit - 500 pts MDM2 GST Tagged protein	01RA
HTRF MDM2 Binding Kit - 500 pts MI-1061 Red Ligand	01RA

Results :

Incubation : 1h 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	4716
Std 1	0,128	4435
Std 2	0,64	4363
Std 3	3,2	3529
Std 4	16	1727
Std 5	80	754
Std 6	400	368
Std 7	2000	387



Assay characteristics :

	Found value	Specifications Rev 01
S/B	13	15-24
IC50	8.2 **	10 - 30 nM

** Batch validated with EC50 a little bit lower than specification

HTRF MDM2 Binding Kit - 500 pts - Lot n° : 01E

Accepted

Rejected

Quality control date :	2024-12-23	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