

HTRF® Research Reagents

HTRF cIAP1 BIR3 Bind Kit - 500 pts

Product Number : 64BDIAPB3PEG

Lot Number : 02B

Expiration date : 2025-09-15

Document version : 1

Reagents :		Lot n°
HTRF Diluent 9 (5X) - 2 mL		04A
HTRF PROTAC Binding Buffer 2 - 20 mL		05RA
HTRF cIAP1 BIR3 Bind Kit - Std		01RA
HTRF cIAP1 BIR3 Bind Kit - 500 pts Tb Cryptate antibody		02RA
HTRF cIAP1 BIR3 Bind Kit - 500 pts cIAP1 BIR3 GST Tagged		02RA
HTRF cIAP1 BIR3 Bind Kit - 500 pts LCL161 Red Ligand		02RA

Results :

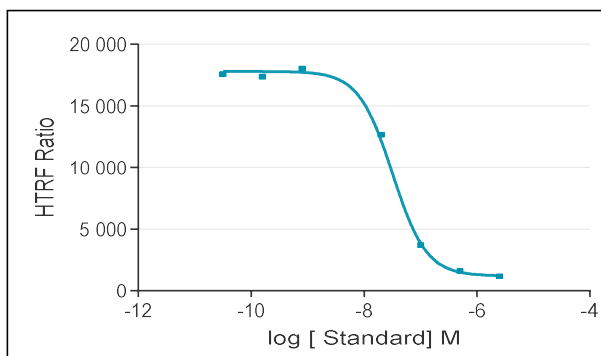
Incubation : 2h at RT

Plate : 384 wells small volume white plate 20 µL

Reader : Pherastar Plus

Fitting model : log(inhibitor) vs.response-Variable slope (four parameters)

cAMP	Conc. (nM)	RATIO
Std 0	0	17590
Std 1	0,16	17394
Std 2	0,8	18029
Std 3	4	20173
Std 4	20	12664
Std 5	100	3719
Std 6	500	1604
Std 7	2500	1176



Assay characteristics :

	Found value	Specifications Rev 01
négative	17590	7129-29023
S/B	15	7-25
IC 50	31,9 nM	12-38 nM

HTRF cIAP1 BIR3 Bind Kit - 500 pts lot 02B

☒ Accepted☐ Rejected

Quality control date : 2024-10-31

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.