

HTRF® Research Reagents

HTRF Human VHL Bind. Kit - 500 pts

Product Number : 64BDVHLPEG Lot Number : 02D

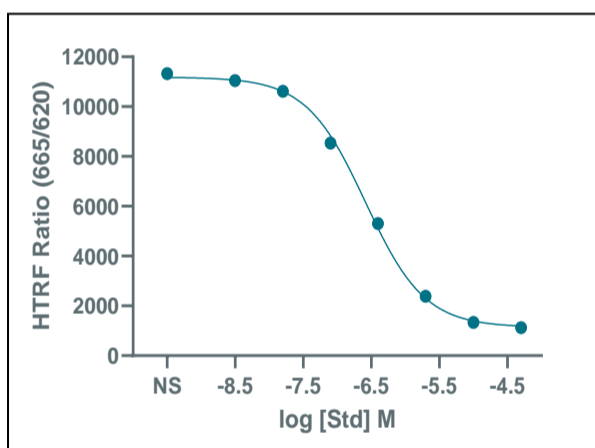
Expiration date : 2025-01-08 Document version : 1

Reagents :	Lot n°
HTRF Diluent 9 (5X) - 2 mL	03B
HTRF PROTAC Binding Buffer 1 - 20 mL	05AR
HTRF Human VHL Bind. Kit - Std	02A
HTRF Human VHL Bind. Kit - 500 pts Eu Cryptate antibody	02RA
HTRF Human VHL Bind. Kit - 500 pts 6His-VHL complex	02RA
HTRF Human VHL Bind. Kit - 500 pts VHL-Red Ligand	02RA

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	11323
Std 1	0,0032	11047
Std 2	0,016	10613
Std 3	0,08	8537
Std 4	0,4	5301
Std 5	2	2389
Std 6	10	1339
Std 7	50	1122



Assay characteristics :

	Found value	Specifications Rev 01
S/B	10	6-12
IC 50	260 nM	150-370 nM

HTRF Human VHL Bind. Kit - 500 pts - Lot n° : 02D

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.