

HTRF® Research Reagents

HTRF Human VHL Bind. Kit - 500 pts

Product Number : 64BDVHLPEG Lot Number : 02E

Expiration date : 2026-01-04 Document version : 1

Reagents :	Lot n°
HTRF Diluent 9 (5X) - 2 mL	03RB
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	03AR
HTRF Human VHL Bind. Kit - 500 pts VHL-Red Ligand	02RA

Results :

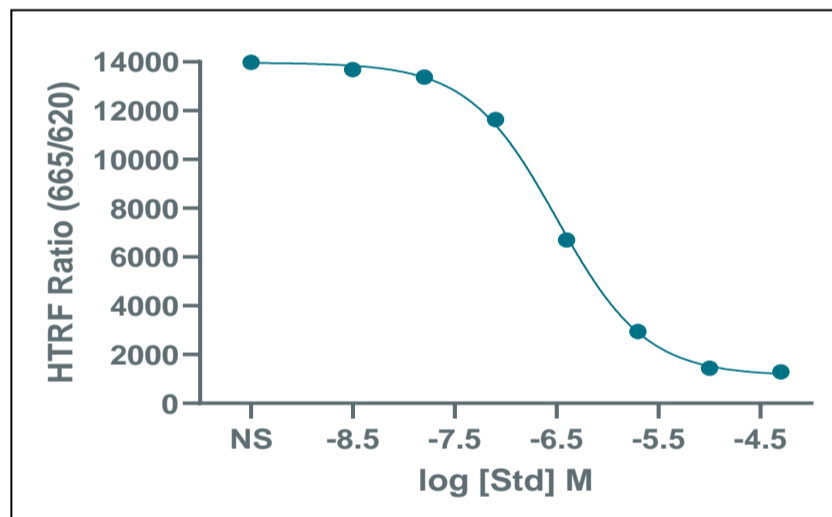
Incubation :

Plate :

Reader : Pherastar Plus

Fitting model :

STD	Conc. (µM)	RATIO
Std 0	0	13983
Std 1	0.0032	13690
Std 2	0.016	13381
Std 3	0.08	11634
Std 4	0.4	6699
Std 5	2	2946
Std 6	10	140
Std 7	50	1296



Assay characteristics :

	Found value	Specifications Rev 01
S/B	11	6-12
IC50	320	150-340

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

Accepted

Rejected

Quality control date : 2024-07-23	Supervisor's initials : ML
-----------------------------------	----------------------------

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