

HTRF® Research Reagents

HTRF Human VHL Bind. Kit - 500 pts

Product Number : 64BDVHLPEG Lot Number : 02F

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		04A
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 :

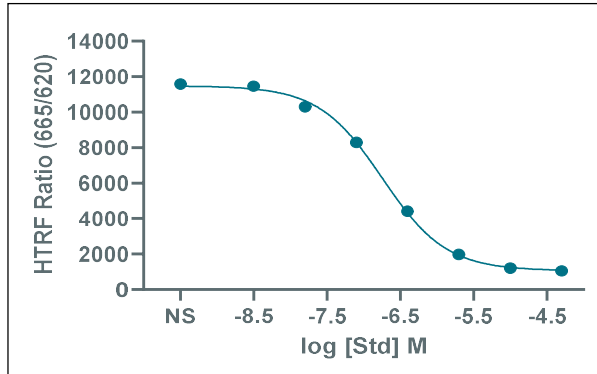
Incubation :

Plate :

Reader : Pherastar Plus

Fitting model :

STD	Conc. (µM)	RATIO
Std 0	0	11585
Std 1	0,0032	11461
Std 2	0,016	10304
Std 3	0,08	8286
Std 4	0,4	4408
Std 5	2	1985
Std 6	10	1212
Std 7	50	1056



Assay characteristics :

	Found value	Specifications Rev 01
S/B	11	6-12
IC50	180	150-340

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

Accepted

Rejected

Quality control date : 2024-09-18	Supervisor's initials : MP
-----------------------------------	----------------------------

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