

HTRF® Research Reagents

HTRF PD1/PD-L1 Bind. kit - 500 pts

Product Number : 64PD1PEG Lot Number : 07D

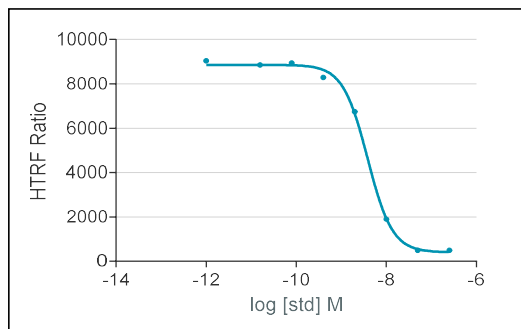
Expiration date : 2026-04-24 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	07RA
HTRF PD1/PD-L1 Bind. kit - 500 pts Tag2-PD1	06RA
HTRF PD1/PD-L1 Bind. kit - Std	05A
HTRF PD1/PD-L1 Bind. kit - 500 pts Eu Cryptate reagent	07A
HTRF PD1/PD-L1 Bind. kit - 500 pts Tag1-PD-L1	06RA
HTRF PD1/PD-L1 Bind. kit - 500 pts XL665 antibody	07A

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

PD1/PD- L1	Conc. (nMf)	RATIO
Std 0	0	9044
Std 1	0,016	8854
Std 2	0,08	8943
Std 3	0,4	8299
Std 4	2	6747
Std 5	10	1901
Std 6	50	494
Std 7	250	499



Assay characteristics :

	Found value	Specifications Rev 01
IC 50	2,87	2 - 7 nM
S/B	18	16-26

HTRF PD1/PD-L1 Bind. kit - 500 pts - Lot n° : 07D

Accepted

Rejected

Quality control date : 2024-08-20	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.