

HTRF® Research Reagents

HTRF S1/ACE2 Bind. kit - 500 pts

Product Number : 64BDS1PEG Lot Number : 02A

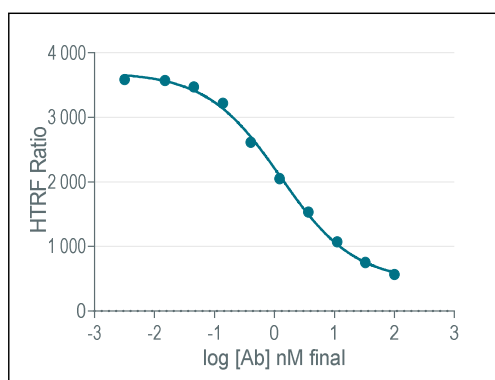
Expiration date : 2025-08-08 Document version : 1

Reagents :		Lot n°
HTRF Detect. Buf. 3 - 50 mL		29RA
HTRF S1/ACE2 Bind. kit - 500 pts d2 reagent		01BR
HTRF S1/ACE2 Bind. kit - 500 pts Eu Cryptate reagent		02AR
HTRF S1/ACE2 Bind. kit - Ctrl		02A
HTRF S1/ACE2 Bind. kit - modop		2

Results :

Incubation : Overnight at 4°C
Plate : 384 wells small volume white plate 20 µL
Reader : Pherastar Plus
Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

Human STING	Conc. (nMi)	RATIO %
Std 0	0	3587
Std 1	0.02	3570
Std 2	0.05	3473
Std 3	0.14	3218
Std 4	0.41	2614
Std 5	1.23	2050
Std 6	3.70	1532
Std 7	11.11	1071
Std 8	33	753
Std 9	100	566



Assay characteristics :

	Found value	Specifications Rev 01
ratio max	3592	3005 - 4510
Negative	339	306 - 460
IC 50	1.2	0.54 - 2 nM

HTRF S1/ACE2 Bind. kit - 500 pts - Lot n° : 02A

☒ Accepted☐ Rejected

Quality control date : 2023-09-27 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.