

HTRF® Research Reagents

HTRF SARS-CoV2 capsid kit - 10K pts

Product Number : 63ADK113PEH Lot Number : 05B

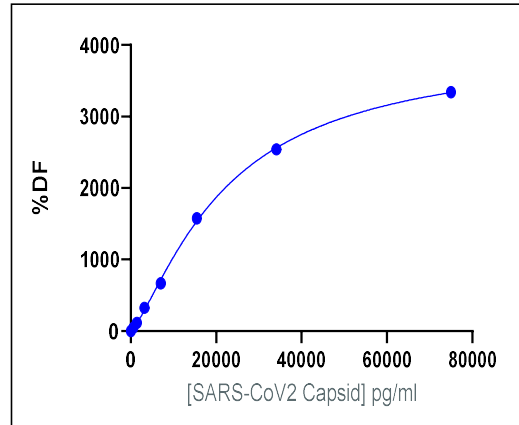
Expiration date : 2025-11-24 Document version : 1

Reagents :	Lot n°
HTRF SARS-CoV2 capsid kit - 10K pts d2 antibody	05RA
HTRF SARS-CoV2 capsid kit - 10K pts Eu Cryptate antibody	05RA
HTRF SARS-CoV2 capsid kit - Std	04RA
HTRF P-T prot. - Lysis Buf.5 (4X) 130 mL	02RA
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

Results :

Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader :

	Conc (pg/ml)	Ratio	RATIO Accepted Range
Std 0	0	538	488 - 535
Std 1	661	855	682 - 870
Std 2	1455	1191	1050 - 1223
Std 3	3202	2117	1850 - 2838
Std 4	7044	4533	4103 - 4656
Std 5	15496	8829	7075 - 11414
Std 6	34091	13548	12879 - 15499
Std 7	75000	17645	16116 - 20060



HTRF SARS-CoV2 capsid kit - 10K pts - Lot n° : 05B

Accepted

Rejected

Quality control date : 2022-12-19	Supervisor's initials : AC
-----------------------------------	----------------------------

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.