

HTRF® Research Reagents

HTRF Viral dsRNA detect. kit - 500 pts

Product Number : 64RNAPEG Lot Number : 09D

Expiration date : 2025-05-05 Document version : 1

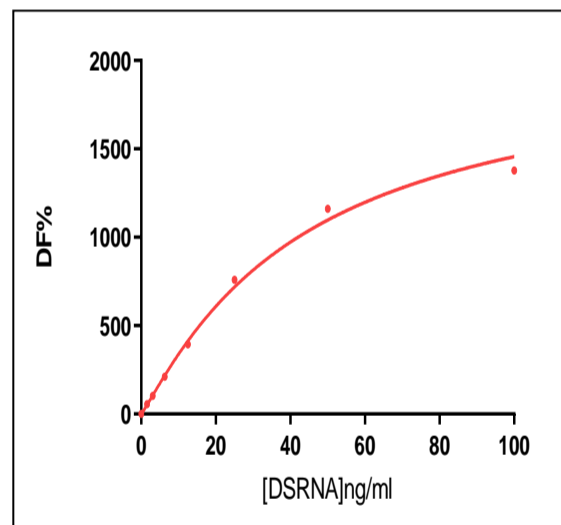
Reagents :	Lot n°
HTRF Detect. Buf. 11 - 7 mL	08AR
HTRF Viral dsRNA detect. kit - 500 pts d2 antibody	09A
HTRF Viral dsRNA detect. kit - 500 pts Eu Cryptate antibody	09A
HTRF Viral dsRNA detect. kit - Std	08RA
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14A

Results :

Incubation : Overnight at 4°C  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : 4 Parameter Logistic (4PL)

Results :

RNA	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0%	0-0
Std 1	1,56	55%	33-62
Std 2	3,13	103%	70-145
Std 3	6,25	211%	145-269
Std 4	12,5	394%	282-523
Std 5	25	759%	492-914
Std 6	50	1161%	748-1389
Std 7	100	1377%	1096-2036



HTRF Viral dsRNA detect. kit - 500 pts - Lot n° : 09D

Accepted

Rejected

Quality control date : 2024-01-11	Supervisor's initials : FMV
-----------------------------------	-----------------------------

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.