

HTRF® Research Reagents

## HTRF Viral dsRNA detect kit - 500 pts

Product Number : 64RNAPEG Lot Number : 08F

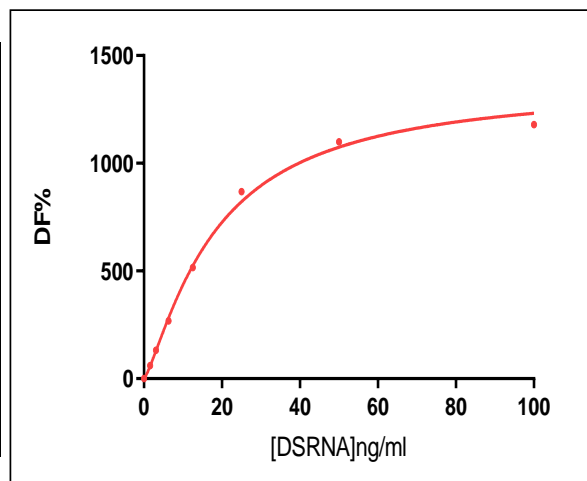
Expiration date : 21/07/2024 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 11 - 7 mL	07RA
HTRF Viral dsRNA detect. kit - 500 pts d2 antibody	08RA
HTRF Viral dsRNA detect. kit - 500 pts Eu Cryptate antibody	08RA
HTRF Viral dsRNA detect. kit - Std	07RA
HTRF Viral dsRNA detect. kit - modop	5
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)

RNA	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0%	0-0
Std 1	1,56	60	33-62
Std 2	3,13	132	70-145
Std 3	6,25	268	145-269
Std 4	12,5	516	282-523
Std 5	25	868	492-914
Std 6	50	1099	748-1389
Std 7	100	1179	1096-2036



HTRF Human Viral DsRNA detect kit - 500 pts - Lot n° : 08F

Accepted

Rejected

Quality control date : 2023-09-13	Supervisor's initials : FMV
-----------------------------------	-----------------------------

The information provided in this document is valid for the specimen 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