

HTRF® Research Reagents

HTRF Viral dsRNA detect. kit - 10K pts

Product Number : 64RNAPEH Lot Number : 06B

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

Reagents :	Lot n°
HTRF Detect. Buf. 11 - 105 mL	06RA
HTRF Viral dsRNA detect. kit - 10K pts d2 antibody	06RA
HTRF Viral dsRNA detect. kit - 10K pts Eu Cryptate antibody	06RA
HTRF Viral dsRNA detect. kit - Std	08RA
HTRF P-T prot. - Lysis Buf.4 (4X) 130 mL	14RA

Results :

Incubation : Overnight at 4°C
 Plate : 384 wells small volume white plate 20 µL
 Reader : /
 Fitting model : /

Assay characteristics :

RNA	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 0	0	0%	0-0
Std 1	1.56	35%	33-62
Std 2	3.13	70%	70-145
Std 3	6.25	151%	145-269
Std 4	12.5	295%	282-523
Std 5	25	602%	492-914
Std 6	50	1113%	748-1389
Std 7	100	1472%	1096-2036

HTRF Viral dsRNA detect. kit - 10K pts - Lot n° : 06B

Accepted

Rejected

Quality control date : 2023-09-22	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.