

HTRF® Research Reagents

HTRF dsRNA (IVT) Kit - 500 pts

Product Number : 64DSRNAPEG Lot Number : 01C

Expiration date : 2028-01-16 Document version : 1

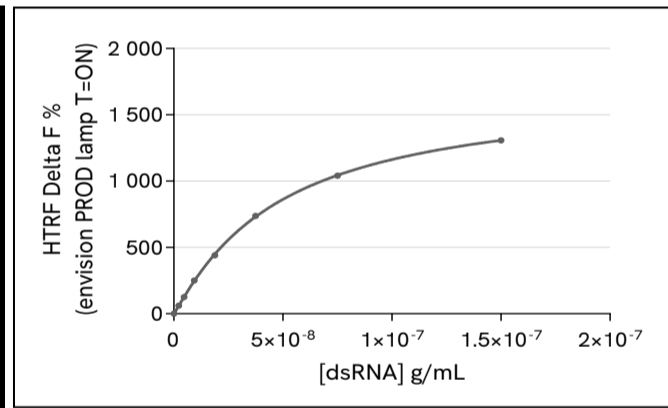
Reagents :	Lot n°
Diluent 14 (4X) - 2 mL	01B
HTRF Detect. Buf. 19 - 7 mL	01A
HTRF dsRNA (IVT) Kit - Std	02A
HTRF dsRNA (IVT) Kit - 500 pts d2 antibody	01B
HTRF dsRNA (IVT) Kit - 500 pts Eu Cryptate antibody	01B

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Envision Lamp Prod
 Fitting model : 4 Parameter Logistic (4PL with 1/Y²)

Assay characteristics :

dsRNA	Conc. (ng/ml)	DF %	DF% Accepted Range
Std 1	2,34	61%	51-86
Std 2	4,69	125%	102-159
Std 3	9,375	251%	215-323
Std 4	18,75	441%	396-594
Std 5	37,5	738%	697-1045
Std 6	75	1041%	1010-1514
Std 7	150	1308%	1247-1847



	Found value	Specifications Rev 01
LoQ	0,8	< 1ng/ml

HTRF dsRNA (IVT) Kit - 500 pts - Lot n° : 01C

Accepted

Rejected

Quality control date : 2025-02-07	Supervisor's initials : PL
-----------------------------------	----------------------------

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.