

## HTRF® Research Reagents

## HTRF dsRNA (IVT) Kit - 500 pts

Product Number : 64DSRNAPEG

Lot Number : 02A

Expiration date : 2028-01-31

Document version :

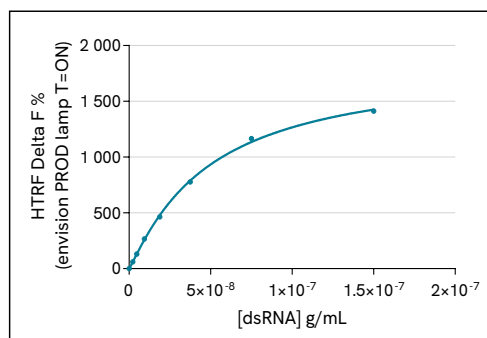
| Reagents :                                          |  | Lot n° |
|-----------------------------------------------------|--|--------|
| Diluent 14 (4X) - 2 mL                              |  | 01B    |
| HTRF Detect. Buf. 19 - 7 mL                         |  | 05A    |
| HTRF dsRNA (IVT) Kit - Std                          |  | 02A    |
| HTRF dsRNA (IVT) Kit - 500 pts d2 antibody          |  | 02A    |
| HTRF dsRNA (IVT) Kit - 500 pts Eu Cryptate antibody |  | 02A    |
|                                                     |  |        |

## 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/Y<sup>2</sup>)

## Assay characteristics :

| dsRNA | Conc.<br>(ng/ml) | DF %  | DF%<br>Accepted<br>Range |
|-------|------------------|-------|--------------------------|
| Std 1 | 2,34             | 61%   | 51-86                    |
| Std 2 | 4,69             | 131%  | 102-159                  |
| Std 3 | 9,375            | 266%  | 215-323                  |
| Std 4 | 18,75            | 464%  | 396-594                  |
| Std 5 | 37,5             | 778%  | 697-1045                 |
| Std 6 | 75               | 1166% | 1010-1514                |
| Std 7 | 150              | 1412% | 1247-1847                |



|                                                                          | Found value | Specifications Rev 01      |
|--------------------------------------------------------------------------|-------------|----------------------------|
| LoD                                                                      | 0,15        | < 1ng/ml                   |
| HTRF dsRNA (IVT) Kit - 500 pts - Lot n° : 02A                            |             |                            |
| <input checked="" type="radio"/> Accepted <input type="radio"/> Rejected |             |                            |
| Quality control date :                                                   | 2025-04-30  | Supervisor's initials : FL |

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.