

## HTRF® Research Reagents

## HTRF Human IL1 beta kit - 96 pts

Product Number : 62HIL1BPET Lot Number : 11B

Expiration date : 2027-08-10

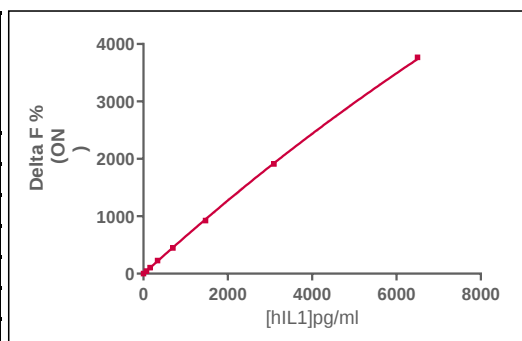
Document version : 1

| Reagents :                                            |  | Lot n° |
|-------------------------------------------------------|--|--------|
| PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND                |  | 23     |
| HTRF Diluent 5 (5X) - 2 mL                            |  | 22A    |
| HTRF Detect. Buf. 3 - 0.5 mL                          |  | 30AR   |
| HTRF Human IL1 beta kit - Std                         |  | 05RA   |
| HTRF Human IL1 beta kit - 96 pts Eu Cryptate antibody |  | 11A    |
| HTRF Human IL1 beta kit - 96 pts XL665 antibody       |  | 11A    |
|                                                       |  |        |

## Results :

Incubation : Overnight at RT  
Plate : 384 wells small volume white plate 20 µL  
Reader : EnVision  
Fitting model : 4 Parameter Logistic (4PL 1/y2)\* model

| IL1 beta | Conc.<br>(pg/ml) | DF %  | DF<br>Accepted<br>Range |
|----------|------------------|-------|-------------------------|
| Std 1    | 75,8             | 46%   | 39-102                  |
| Std 2    | 159,1            | 106%  | 80-219                  |
| Std 3    | 334,2            | 228%  | 175-403                 |
| Std 4    | 701,9            | 447%  | 382-793                 |
| Std 5    | 1474             | 926%  | 837-1508                |
| Std 6    | 3095             | 1910% | 1704-2800               |
| Std 7    | 6500             | 3769% | 2840-4975               |



## Assay characteristics :

|                  | Found value | Specifications Rev 02 |
|------------------|-------------|-----------------------|
| Detection limite | 31 pg/ml    | ≤ 75.8 pg/ml          |

HTRF Human IL1 beta kit - 96 pts - Lot n° : 11B

☒ Accepted☐ Rejected

|                        |            |                         |    |
|------------------------|------------|-------------------------|----|
| Quality control date : | 2025-04-28 | 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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.