



HTRF® Research Reagents

HTRF Human IL6 kit - 96 pts

Product Number :62HIL06PET

Lot Number :12B

Expiration date :2027-02-07

Document version :1

| Reagents :                                       | Lot n° |
|--------------------------------------------------|--------|
| PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND           | 2413   |
| HTRF Diluent 5 (5X) - 2 mL                       | 21B    |
| HTRF Detect. Buf. 3 - 0.5 mL                     | 30AR   |
| HTRF Human IL6 kit - Std                         | 06RA   |
| HTRF Human IL6 kit - 96 pts Eu Cryptate antibody | 11A    |
| HTRF Human IL6 kit - 96 pts XL665 antibody       | 12A    |
|                                                  |        |

Results :

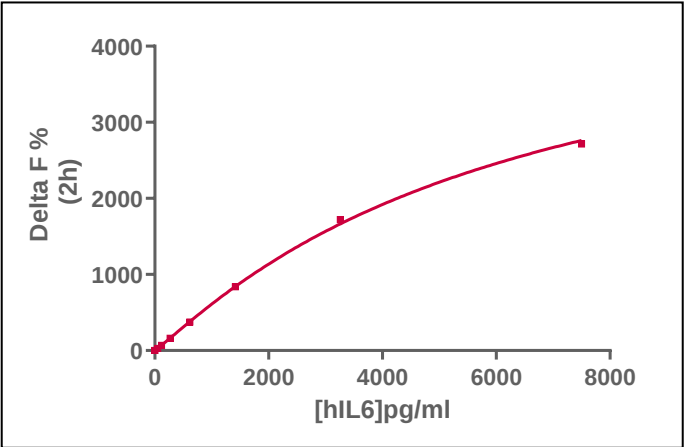
Incubation : 2h at RT

Plate : 384 wells small volume white plate 20 µL

Reader : Envision

Fitting model : 4 Parameter Logistic (4PL 1/y2)\* model

| hIL6  | Conc.<br>(pg/ml) | DF %  | DF<br>Accepted<br>Range |
|-------|------------------|-------|-------------------------|
| Std 1 | 50,7             | 28%   | 25-57                   |
| Std 2 | 117              | 67%   | 60-118                  |
| Std 3 | 268              | 163%  | 128-251                 |
| Std 4 | 616              | 374%  | 300-495                 |
| Std 5 | 1417             | 837%  | 632-1034                |
| Std 6 | 3260             | 1723% | 1255-2095               |
| Std 7 | 7500             | 2714% | 1875-3131               |



Assay characteristics :

|                 | Found value | Specifications Rev 03 |
|-----------------|-------------|-----------------------|
| Detection limit | 7,3 pg/ml   | ≤ 50.7 pg/ml          |

HTRF Human IL6 kit - 96 pts - Lot n° : 12B

Accepted

Rejected

|                                  |                           |
|----------------------------------|---------------------------|
| Quality control date :2025-01-22 | 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.

