

HTRF® Research Reagents

HTRF Human IL17 kit - 500 pts

Product Number : 62HIL17PEG Lot Number : 04B

Expiration date : 2024/07/21 Document version : 1

| Reagents :   | Lot n° |
|--|--------|
| HTRF Diluent 5 (5X) - 2 mL                         | 21RA   |
| HTRF Detect. Buf. 3 - 1.5 mL                       | 29A    |
| HTRF Human IL17 kit - Std                          | 03RA   |
| HTRF Human IL17 kit - 500 pts d2 antibody          | 04A    |
| HTRF Human IL17 kit - 500 pts Eu Cryptate antibody | 04A    |
| HTRF Human IL17 kit - modop                        | 7      |

Results :

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

| IL17  | Conc. (pg/ml) | DF %  | DF Accepted Range |
|-------|---------------|-------|-------------------|
| Std 1 | 88            | 60%   | 48-66             |
| Std 2 | 194           | 122%  | 99-134            |
| Std 3 | 427           | 245%  | 185-275           |
| Std 4 | 939           | 511%  | 366-612           |
| Std 5 | 2066          | 985%  | 739-1275          |
| Std 6 | 4545          | 1885% | 1556-2551         |
| Std 7 | 10000         | 3136% | 2676-3186         |

Assay characteristics :

|                 | Found value | Specifications Rev 02 |
|-----------------|-------------|-----------------------|
| Detection limit | 6 pg/ml     | ≤ 8 pg/ml             |
|                 | /           | /                     |
|                 | /           | /                     |
|                 | /           | /                     |

HTRF Human IL17 kit - 500 pts - Lot n° : 04B

Accepted

Rejected

|                                   |                            |
|-----------------------------------|----------------------------|
| Quality control date : 2023-09-19 | 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.