

HTRF® Research Reagents

HTRF (h) IL12/23 p40 Kit - 500 pts

Product Number : 62HIL12P40PEG Lot Number : 01A

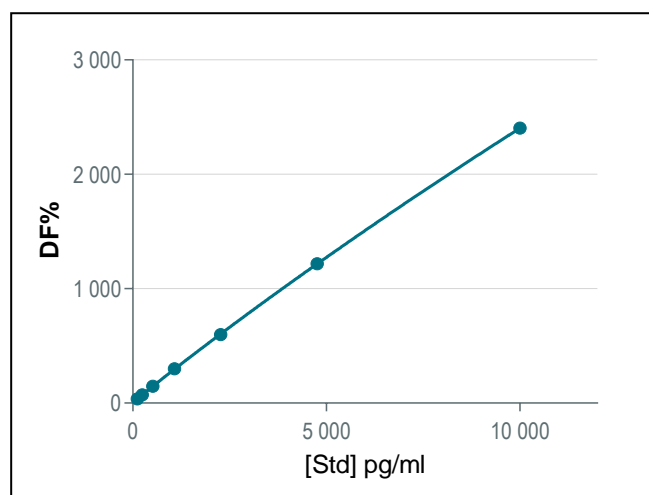
Expiration date : 2025-10-28 Document version : 1

| Reagents : | Lot n° |
|---|--------|
| HTRF Diluent 5 (5X) - 2 mL | 21A |
| HTRF Detect. Buf. 3 - 1.5 mL | 28A |
| HTRF (h) IL12/23 p40 Kit - Std | 01A |
| HTRF (h) IL12/23 p40 Kit - 500 pts d2 antibody | 01A |
| HTRF (h) IL12/23 p40 Kit - 500 pts Eu Cryptate antibody | 01A |
| HTRF (h) IL12/23 p40 Kit - modop | 1 |

Results :

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

| h IL12P40 Kit | Conc. (pg/ml) | DF % | DF Accepted Range |
|---------------|---------------|------|-------------------|
| Std 1 | 117 | 37 | 28 --- 40 |
| Std 2 | 245 | 72 | 52 --- 80 |
| Std 3 | 514 | 147 | 118 --- 159 |
| Std 4 | 1080 | 300 | 235 --- 310 |
| Std 5 | 2268 | 598 | 541 --- 607 |
| Std 6 | 4762 | 1218 | 1078 --- 1262 |
| Std 7 | 10000 | 2404 | 2162 --- 2528 |



Assay characteristics :

| | Found value | Specifications Rev 01 |
|-----------------|-------------|-----------------------|
| Detection limit | 12.6 pg/ml | < 117 pg/ml |

HTRF (h) IL12/23 p40 Kit - 500 pts - Lot n° : 01A

Accepted

Rejected

| | |
|-----------------------------------|-----------------------------|
| Quality control date : 2024-07-09 | Supervisor's initials : FMV |
|-----------------------------------|-----------------------------|

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.