

HTRF® Research Reagents

HTRF Human IFN g kit - 96 pts

Product Number : 62HIFNGPET Lot Number : 10L

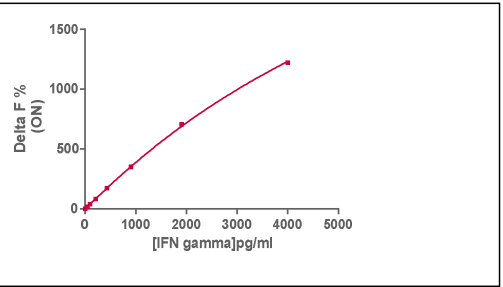
Expiration date : 2024/12/14 Document version : 1

| Reagents : | Lot n° |
|--|--------|
| PLAQUE BLANCHE 96 PUIFS SV COSMOBRAND | 2203 |
| HTRF Diluent 5 (5X) - 2 mL | 21RA |
| HTRF Detect. Buf. 3 - 0.5 mL | 28RA |
| HTRF Human IFN g kit - Std | 11RA |
| HTRF Human IFN gamma kit - modop | 9 |
| HTRF Human IFN g kit - 96 pts Eu Cryptate antibody | 10A |
| HTRF Human IFN g kit - 96 pts XL665 antibody | 10A |

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**

| hIFNg | Conc. (pg/ml) | DF % | DF Accepted Range |
|-------|---------------|-------|-------------------|
| Std 1 | 46.6 | 18% | 13-21 |
| Std 2 | 98 | 39% | 25-48 |
| Std 3 | 205.7 | 81% | 60-96 |
| Std 4 | 432 | 172% | 129-195 |
| Std 5 | 907 | 351% | 263-387 |
| Std 6 | 1904.8 | 704% | 556-747 |
| Std 7 | 4000 | 1220% | 1026-1310 |



Assay characteristics :

| | Found value | Specifications Rev 01 |
|--|------------------------|-----------------------|
| | Detection limit | 10 pg/ml |
| | / | / |
| | / | / |
| | / | / |

HTRF Human IFN g kit - 96 pts - Lot n° : 10L

Accepted

Rejected

| | |
|-----------------------------------|----------------------------|
| Quality control date : 2023-09-18 | 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.