

HTRF® Research Reagents

### HTRF Human IFN g kit - 96 pts

Product Number : 62HIFNGPET Lot Number : 11B

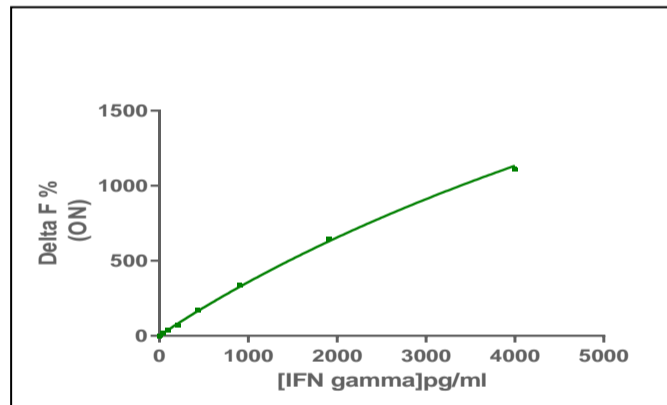
Expiration date : 2027-01-09 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 0.5 mL	30AR
HTRF Human IFN g kit - Std	12A
HTRF Human IFN g kit - 96 pts Eu Cryptate antibody	08RB
HTRF Human IFN g kit - 96 pts XL665 antibody	11RB

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	20%	13-21
Std 2	98	40%	25-48
Std 3	205,7	71%	60-96
Std 4	432	171%	129-195
Std 5	907	336%	263-387
Std 6	1904,8	646%	556-747
Std 7	4000	1110%	1026-1310



Assay characteristics :

	Found value	Specifications Rev 01
<b>Detection limit</b>	<b>9,5 pg/ml</b>	<b>≤ 46.6 pg/ml</b>

HTRF Human IFN g kit - 96 pts - Lot n° : 11B

Accepted

Rejected

Quality control date : 2024-03-27	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.