

HTRF® Research Reagents

HTRF Human IFN g kit - 500 pts

Product Number : 62HIFNGPEG Lot Number : 12J

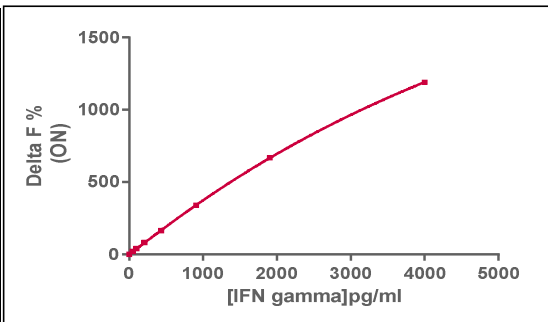
Expiration date : 2025-08-17 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	28A
HTRF Human IFN g kit - Std	11RA
HTRF Human IFN g kit - 500 pts Eu Cryptate antibody	11RB
HTRF Human IFN g kit - 500 pts XL665 antibody	12RA
HTRF Human IFN gamma kit - modop	8

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	21%	13-21
Std 2	98	40%	25-48
Std 3	205.7	82%	60-96
Std 4	432	164%	129-195
Std 5	907	340%	263-387
Std 6	1904.8	668%	556-747
Std 7	4000	1191%	1026-1310



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	9 pg/ml	< 46.6 pg/ml

HTRF Human IFN g kit - 500 pts - Lot n° : 12J

Accepted

Rejected

Quality control date : 2023-11-17	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.