

HTRF® Research Reagents

HTRF Human IFN g kit - 10K pts

Product Number : 62HIFNGPEH Lot Number : 13A

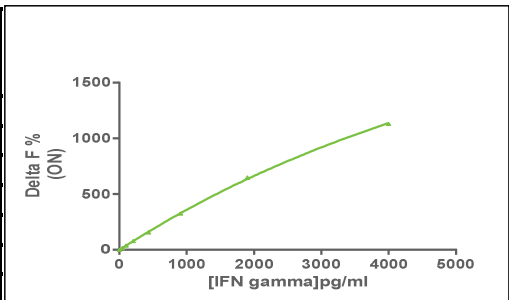
Expiration date : 2025-08-17 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Human IFN g kit - 10K pts Eu Cryptate antibody	13AR
HTRF Human IFN g kit - 10K pts XL665 antibody	13AR
HTRF Human IFN g kit - Std	11RA
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	17%	13-21
Std 2	98	38%	25-48
Std 3	205.7	78%	60-96
Std 4	432	158%	129-195
Std 5	907	325%	263-387
Std 6	1904.8	649%	556-747
Std 7	4000	1130%	1026-1310



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	11 pg/ml	< 46.6 pg/ml
	/	/
	/	/
	/	/

HTRF Human IFN g kit - 10K pts - Lot n° : 13A

Accepted

Rejected

Quality control date : 2023-10-12	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.