

HTRF® Research Reagents

HTRF Human IFN alpha kit - 500 pts

Product Number : 62HIFNAPEG Lot Number : 05J

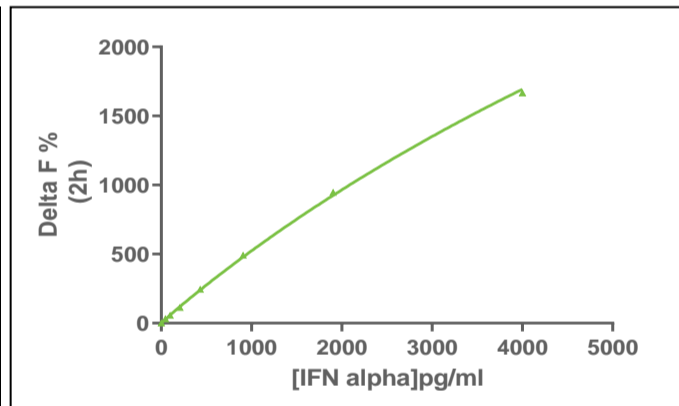
Expiration date : 2025-03-20 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Human IFN alpha kit - Std	05RA
HTRF Human IFN alpha kit - 500 pts d2 antibody	05RA
HTRF Human IFN alpha kit - 500 pts Eu Cryptate antibody	05B

Results :

Incubation : **2h at RT**
 Plate : **384 wells small volume white plate 20 µL**
 Reader : **Pherastar Plus**
 Fitting model : **4 Parameter Logistic (4PL 1/y2)* model**

hIFN alpha	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	47	30%	22-32
Std 2	98	55%	43-66
Std 3	206	113%	106-122
Std 4	432	244%	199-277
Std 5	907	490%	386-537
Std 6	1905	945%	765-961
Std 7	4000	1666%	1247-1551



*This batch has been validated despite std7 out of range

Assay characteristics :

	Found value	Specifications Rev 01
Detection Limit	6 pg/ml	< 47 pg/ml

HTRF Human IFN alpha kit - 500 pts - Lot n° : 05J

Accepted

Rejected

Quality control date : 2024-07-01	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.