

Quality Control report

Human IFN alpha kit - 500 pts
(62HIFNAPEG)

Lot n°06D

Expiration date : 2024-12-21

Reagents :	Lot n°
Diluent 5 (5X) - 2 ml	19G
Human IFN alpha kit - modop	7
Human IFN alpha kit - Standard	05A
Human IFN alpha kit - 500 pts d2 antibody	05A
Human IFN alpha kit - 500 pts Eu Cryptate antibody	06A
Detection buffer 3 - 1.5 mL	28A

Results :

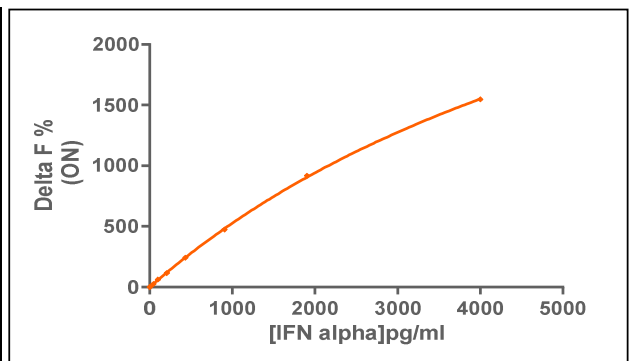
Incubation : 2h at RT

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	26%	22-32
Std 2	98	63%	43-66
Std 3	206	116%	106-122
Std 4	432	242%	199-277
Std 5	907	474%	386-537
Std 6	1905	918%	765-961
Std 7	4000	1548%	1247-1551



Assay characteristics :

	Found value	Specifications
Detection limit	7.2 pg/ml	< 47 pg/ml

Human IFN alpha kit - 500 pts Lot n°06D

Accepted

Rejected

Quality control date :	2022-09-26	Supervisor's initials :	PL
------------------------	------------	-------------------------	----

