

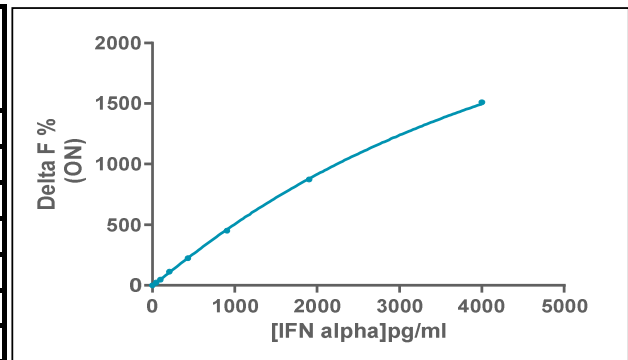
Quality Control report
 Human IFN alpha kit - 500 pts
 (62HIFNAPEG)
 Lot n°06E
 Expiration date : 2024-12-21

Reagents :	Lot n°
Diluent 5 (5X) - 2 ml	21A
Human IFN alpha kit - modop	7
Human IFN alpha kit - Standard	04A
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	22%	22-32
Std 2	98	48%	43-66
Std 3	206	112%	106-122
Std 4	432	224%	199-277
Std 5	907	450%	386-537
Std 6	1905	874%	765-961
Std 7	4000	1510%	1247-1551



Assay characteristics :

	Found value	Specifications
Detection limit	8.9 pg/ml	< 47 pg/ml

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

Accepted

Rejected