



Quality Control report

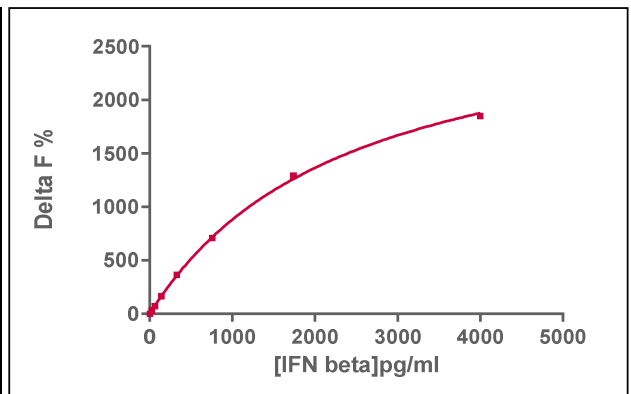
HTRF Human IFN beta kit - 500 pts
 (62HIFNBPEG)
 Lot n°05C
 Expiration date : 2025-11-05

Reagents :	Lot n°
HTRF Activation reagent - 1 mL	03A
HTRF Diluent 5 (5X) - 2 mL	21A
HTRF Detect. Buf. 3 - 1.5 mL	28A
HTRF Human IFN beta kit - Std	05A
HTRF Human IFN beta kit - 500 pts d2 antibody	05RA
HTRF Human IFN beta kit - 500 pts Eu Cryptate antibody	05RA
HTRF Human IFN beta kit - modop	4

Results :

Incubation : 3h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL 1/y2)* model

hIFN beta	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	27	40%	19--51%
Std 2	62	88%	54--115%
Std 3	143	188%	138--231%
Std 4	329	427%	273--493%
Std 5	756	821%	633--929%
Std 6	1739	1457%	1210--1579%
Std 7	4000	2065%	1773--2085%



Assay characteristics :

	Found value	Specifications
Detection Limite	3 pg/ ml	< 5.3pg/ml

HTRF Human IFN beta kit - 500 pts Lot n°05C

Accepted

Rejected

Quality control date :	2023-07-28	Supervisor's initials :	PL
------------------------	------------	-------------------------	----