



Quality Control report

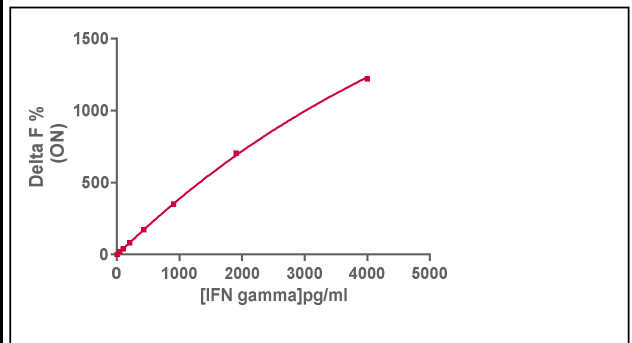
HTRF Human IFN g kit - 96 pts
 (62HIFNGPET)
 Lot n°10G
 Expiration date : 2024-12-14

Reagents :	Lot n°
HTRF Human IFN g kit - Std	11A
HTRF Diluent 5 (5X) - 2 mL	19G
HTRF Detect. Buf. 3 - 0.5 mL	28A
HTRF Human IFN g kit - 96 pts Eu Cryptate antibody	10A
HTRF Human IFN g kit - 96 pts XL665 antibody	10A
HTRF Human IFN g kit - modop	9
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203

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	18%	13-21
Std 2	98	39%	25-48
Std 3	205.7	81%	60-96
Std 4	432	172%	129-195
Std 5	907	351%	263-387
Std 6	1904.8	704%	556-747
Std 7	4000	1220%	1026-1310



Assay characteristics :

	Found value	Specifications
Detection limit	10 pg/ml	≤ 46.6 pg/ml

HTRF Human IFN g kit - 96 pts Lot n°10G

Accepted

Rejected

Quality control date :	2023-08-09	Supervisor's initials :	MP
------------------------	------------	-------------------------	----

