



# Quality Control report

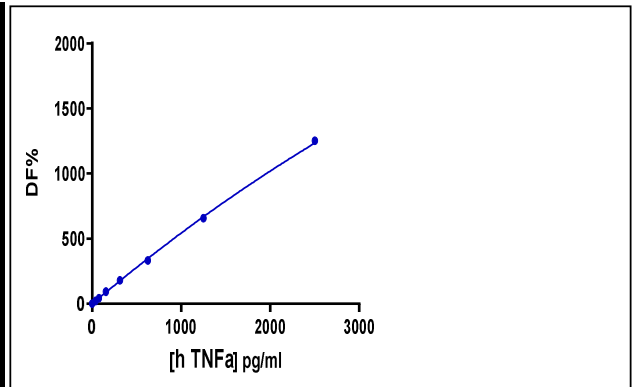
HTRF Human TNF alpha kit - 500 pts  
 ( 62HTNFAPEG )  
 Lot n°09C  
 Expiration date : 2025-09-07

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21A
HTRF Detect. Buf. 3 - 1.5 mL	28A
HTRF Human TNF alpha kit - Std	05A
HTRF Human TNF alpha kit - 500 pts d2 antibody	09RA
HTRF Human TNF alpha kit - 500 pts Eu Cryptate antibody	09RA
HTRF Human TNF alpha kit - modp	7

**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

hTNF alpha	Conc. (pg/ml)	DF %	DF Accepted Range Revision 2
Std 1	39	21%	13-33
Std 2	78	43%	34-61
Std 3	156	92%	70-119
Std 4	312	179%	120-216
Std 5	625	332%	220-359
Std 6	1250	657%	400-645
Std 7	2500	1252%	730-1100



This batch has been validated despite std 6 and 7 out of range

**Assay characteristics :**

	Found value	Specifications
Detection limit	7 pg/ml	< 39 pg/ml

HTRF Human TNF alpha kit - 500 pts Lot n°09C

Accepted

Rejected

Quality control date :	2023-07-20	Supervisor's initials :	PL
------------------------	------------	-------------------------	----

