

Quality Control report

HTRF Human IL2 kit - 96 pts

(62HIL02PET)

Lot n°01X

Expiration date : 2024-03-20

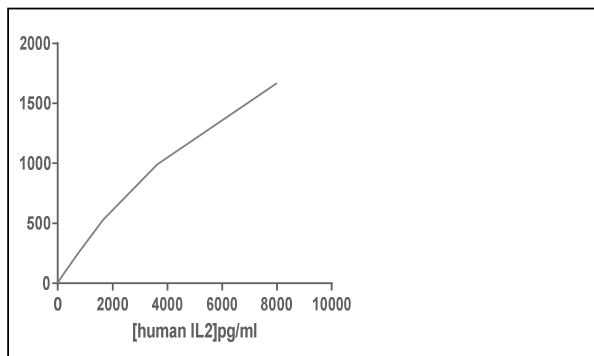
Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 0.5 mL	28A
HTRF Human IL2 kit - Std	07A
HTRF Human IL2 kit - modop	8
HTRF Human IL2 kit - 96 pts d2 antibody	01A
HTRF Human IL2 kit - 96 pts Eu Cryptate antibody	01A

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

Assay characteristics :

IL2	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	70.56	27%	10-45
Std 2	155.23	58%	35-89
Std 3	341.51	121%	100-172
Std 4	751.3	255%	182-363
Std 5	1652.9	530%	389-695
Std 6	3636.4	993%	742-1171
Std 7	8000	1670%	1172-1670



	Found value	Specifications Rev
Detection limit	11 pg/ml	≤ 20 pg/ml

HTRF Human IL2 kit - 96 pts Lot n°01X

Accepted

Rejected

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

