

**Quality Control report**

HTRF HIV p24 kit - 10K pts

( 64P24PEH )

Lot n°03A

Expiration date : 2025-01-22

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	19 A
HTRF Detect. Buf. 3 - 50 mL	28 A
HTRF HIV p24 kit - 10K pts d2 antibody	03A
HTRF HIV p24 kit - 10K pts Eu Cryptate antibody	03A
HTRF HIV p24 kit - Std	02A
HTRF HIV p24 kit - modop	1

**Results :**

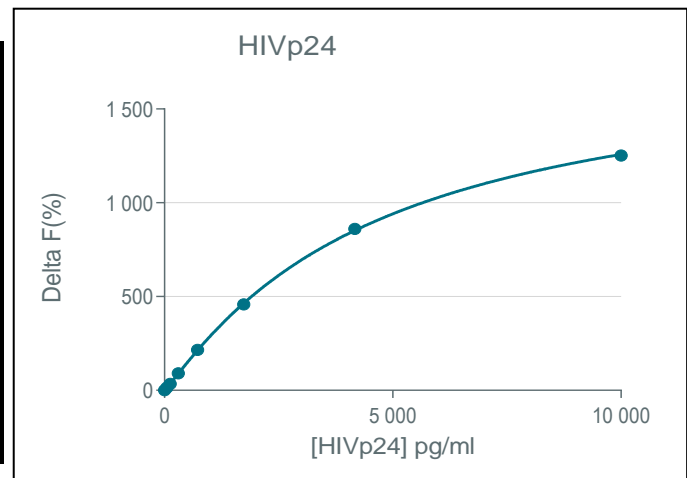
Incubation : 2h at RT

Plate : 384 wells small volume white plate 20 µL

Reader : Pherastar Plus

Fitting model : 4 Parameter Logistic (4PL)

HIV p24	Conc (ng/ml)	Delta F %	Delta F% Accepted Range
Std 1	52	14%	14-21
Std 2	126	34%	28-41
Std 3	301	90%	61-92
Std 4	723	215%	145-217
Std 5	1736	458%	307-460
Std 6	4167	861%	600-900
Std 7	10000	1252%	971-1456



HTRF HIV p24 kit - 10K pts Lot n°03A

Accepted

Rejected

Quality control date :	2022-06-22	Supervisor's initials :	LR
------------------------	------------	-------------------------	----

