

**Quality Control report**

HTRF HIV p24 kit - 10K pts

( 64P24PEH )

Lot n°07B

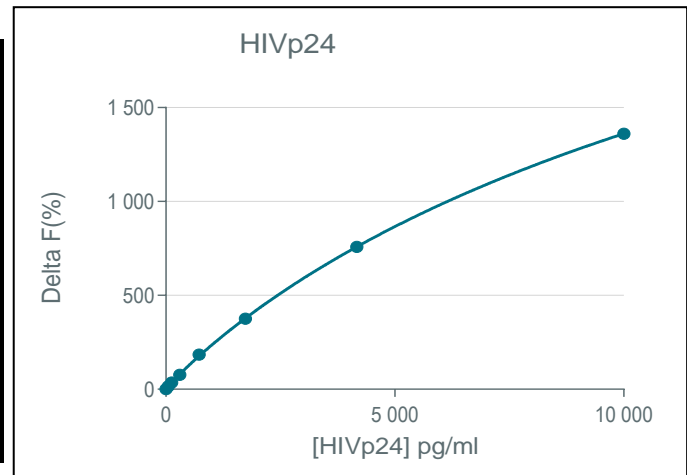
Expiration date : 2025-08-25

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21A
HTRF Detect. Buf. 3 - 50 mL	29A
HTRF HIV p24 kit - 10K pts d2 antibody	05B
HTRF HIV p24 kit - 10K pts Eu Cryptate antibody	07A
HTRF HIV p24 kit - Std	03A
HTRF HIV p24 kit - modop	2

**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	16%	14-21
Std 2	126	35%	28-41
Std 3	301	77%	61-92
Std 4	723	184%	145-217
Std 5	1736	375%	307-460
Std 6	4167	758%	600-900
Std 7	10000	1360%	971-1456



HTRF HIV p24 kit - 10K pts Lot n°07B

Accepted

Rejected

Quality control date :	2023-04-14	Supervisor's initials :	FMV
------------------------	------------	-------------------------	-----

