

Quality Control report

HTRF HIV p24 kit - 10K pts
(64P24PEH)

Lot n°08B

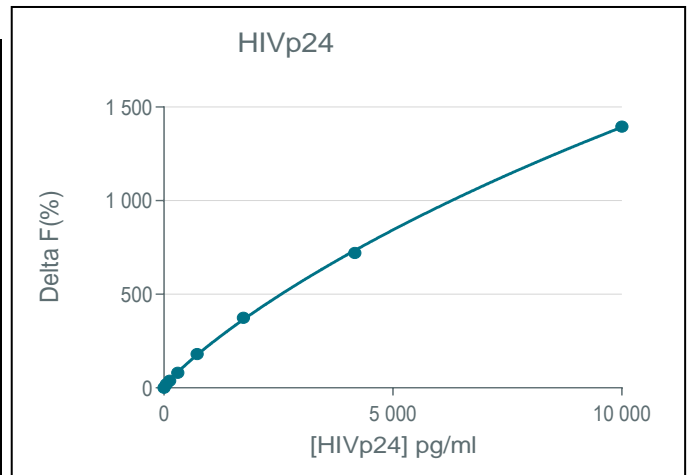
Expiration date : 2026-02-07

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	08A
HTRF HIV p24 kit - 10K pts Eu Cryptate antibody	07A
HTRF HIV p24 kit - Std	04B
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	19%	14-21
Std 2	126	38%	28-41
Std 3	301	81%	61-92
Std 4	723	180%	145-217
Std 5	1736	374%	307-460
Std 6	4167	721%	600-900
Std 7	10000	1395%	971-1456



HTRF HIV p24 kit - 10K pts Lot n°08B

Accepted

Rejected