

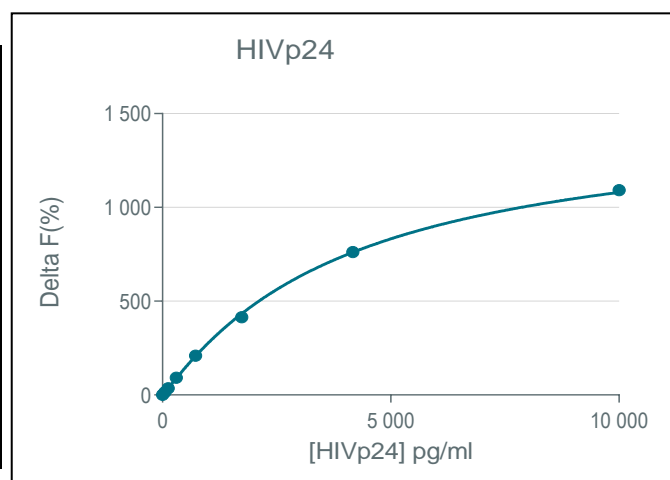
Quality Control report
 HTRF HIV p24 kit - 10K pts
 (64P24PEH)
 Lot n°04A
 Expiration date : 2025-01-22

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	04A
HTRF HIV p24 kit - 10K pts Eu Cryptate antibody	04A
HTRF HIV p24 kit - Std	02A
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	14%	14-21
Std 2	126	36%	28-41
Std 3	301	92%	61-92
Std 4	723	209%	145-217
Std 5	1736	414%	307-460
Std 6	4167	762%	600-900
Std 7	10000	1092%	971-1456



HTRF HIV p24 kit - 10K pts Lot n°04A

Accepted

Rejected

Quality control date : 2022-07-19	Supervisor's initials : FL
-----------------------------------	----------------------------

