

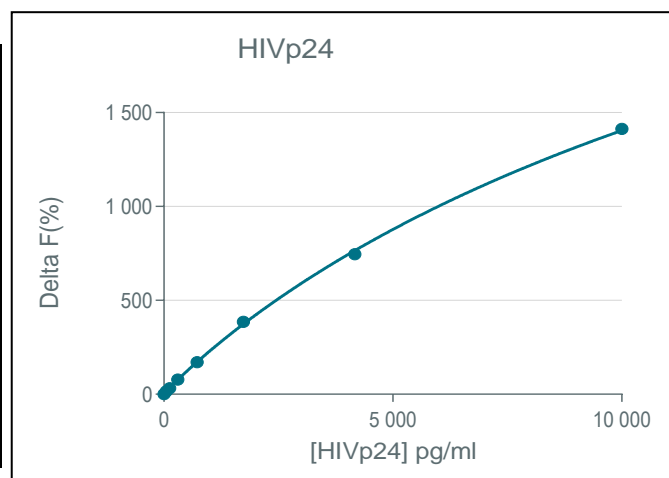
Quality Control report
HTRF HIV p24 kit - 10K pts
(64P24PEH)
Lot n°06A
Expiration date : 2025-01-22

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	19 A
HTRF Detect. Buf. 3 - 50 mL	29A
HTRF HIV p24 kit - 10K pts d2 antibody	06A
HTRF HIV p24 kit - 10K pts Eu Cryptate antibody	06A
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	15%	14-21
Std 2	126	32%	28-41
Std 3	301	78%	61-92
Std 4	723	170%	145-217
Std 5	1736	385%	307-460
Std 6	4167	746%	600-900
Std 7	10000	1413%	971-1456



HTRF HIV p24 kit - 10K pts Lot n°06A

Accepted

Rejected

Quality control date : 2022-12-05	Supervisor's initials : FMV
-----------------------------------	-----------------------------

