



Quality Control report

HTRF HIV p24 kit - 10K pts

(64P24PEH)

Lot n°08A

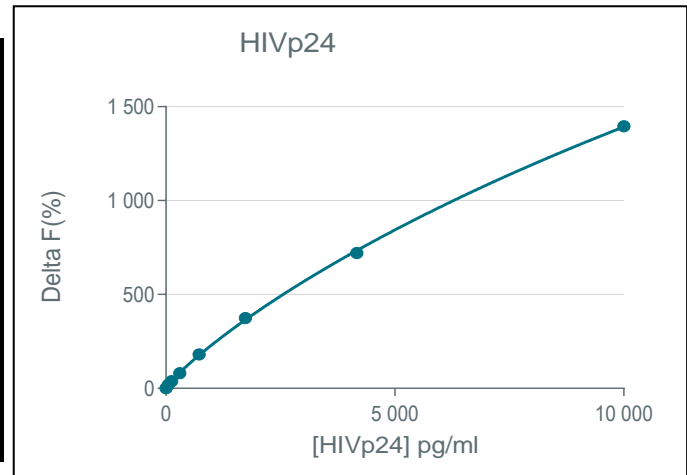
Expiration date : 2026-02-07

| 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°08A

Accepted

Rejected

| | | | |
|------------------------|------------|-------------------------|-----|
| Quality control date : | 2023-06-16 | Supervisor's initials : | FMV |
|------------------------|------------|-------------------------|-----|