

**Quality Control report**

HTRF HIV p24 kit - 10K pts

( 64P24PEH )

Lot n°05A

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          | 05B    |
| HTRF HIV p24 kit - 10K pts Eu Cryptate antibody | 05B    |
| HTRF HIV p24 kit - Std                          | 03A    |
| HTRF HIV p24 kit - modop                        | 2      |

**Results :**

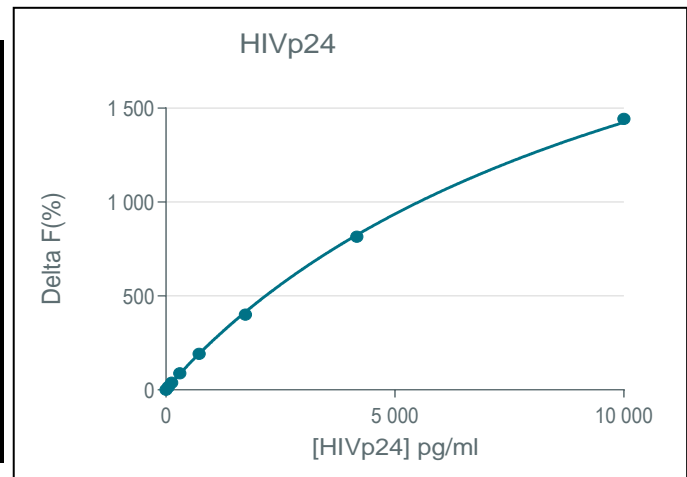
Incubation : 2h at RT

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          | 38%       | 28-41                   |
| Std 3   | 301          | 88%       | 61-92                   |
| Std 4   | 723          | 192%      | 145-217                 |
| Std 5   | 1736         | 400%      | 307-460                 |
| Std 6   | 4167         | 816%      | 600-900                 |
| Std 7   | 10000        | 1443%     | 971-1456                |



HTRF HIV p24 kit - 10K pts Lot n°05A

Accepted

Rejected

|                        |            |                         |     |
|------------------------|------------|-------------------------|-----|
| Quality control date : | 2022-09-29 | Supervisor's initials : | FMV |
|------------------------|------------|-------------------------|-----|

