

ALDO-US

Protocole/Protocol : 3h

Lot N°/Batch # : 317U

| Réactifs Reagents | Lot N° Batch # | Péremption Expiry Date |
|----------------------|-------------------|---------------------------|
| CT-ALDO-RIACT | 142A | 02/03/2022 |
| CAL0-ALDO-RIACT | 26A | 27/04/2023 |
| CAL1-ALDO-RIACT | 25A | 18/11/2022 |
| CAL2-ALDO-RIACT | 25A | 18/11/2022 |
| CAL3-ALDO-RIACT | 25A | 18/11/2022 |
| CAL4-ALDO-RIACT | 25A | 18/11/2022 |
| CAL5-ALDO-RIACT | 25A | 18/11/2022 |
| CONT-ALDO-RIACT | 25A | 18/11/2022 |
| TRA-ALDO-RIACT | 317A | 28/01/2022 |
| ALDO-RIACT MO US | 018 | |

RESULTATS / RESULTS

| Paramètre Parameter | Conc. Conc. | Unité Unit | Réponse Response | Valeur mesurée Measured value |
|------------------------|----------------|---------------|---------------------|----------------------------------|
| CAL.0 | 0.00 | pg/mL | B0/T % | 30.82 |
| CAL.1 | 24.00 | pg/mL | B/B0 | 87.11 |
| CAL.2 | 60.00 | pg/mL | B/B0 | 73.56 |
| CAL.3 | 180.00 | pg/mL | B/B0 | 55.01 |
| CAL.4 | 550.00 | pg/mL | B/B0 | 34.65 |
| CAL.5 | 1 450.00 | pg/mL | B/B0 | 21.34 |

Activité totale / Total activity : 36136 cpm

| Contrôles Controls | Unité Unit | Valeurs attendues Expected values | Valeur mesurée Measured value |
|-----------------------|---------------|--------------------------------------|----------------------------------|
| CONTROL | pg/mL | 70.000 - 130.000 | 81.34 |

Date du contrôle
Control Date

10/11/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

28/01/2022

