

ALDO-US

Protocole/Protocol : 3h

Lot N°/Batch # : 318U

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ALDO-RIACT	318A	24/03/2022
CT-ALDO-RIACT	143A	15/06/2022
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
ALDO-RIACT US MO	018	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	pg/mL	B0/T %	31.82
CAL.1	24.00	pg/mL	B/B0	86.15
CAL.2	60.00	pg/mL	B/B0	72.68
CAL.3	180.00	pg/mL	B/B0	53.22
CAL.4	550.00	pg/mL	B/B0	33.57
CAL.5	1 450.00	pg/mL	B/B0	22.13

Activité totale / Total activity : 38319 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL-	pg/mL	70.000 - 130.000	94.17

Date du contrôle
Control Date

06/01/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

24/03/2022

