

TRYPSS-US

Protocole/Protocol : 3h + 30mn

Lot N°/Batch # : 370U

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCFE07-TRYPSS	370A	21/03/2022
Ab-OCFE07-TRYPSS	24A	07/12/2023
RIP-OCFE07-TRYPSS	20A	25/04/2023
CAL0-OCFE07-TRYPSS	31A	06/11/2023
CAL1-OCFE07-TRYPSS	31A	06/11/2023
CAL2-OCFE07-TRYPSS	31A	06/11/2023
CAL3-OCFE07-TRYPSS	31A	06/11/2023
CAL4-OCFE07-TRYPSS	31A	06/11/2023
CAL5-OCFE07-TRYPSS	31A	06/11/2023
CAL6-OCFE07-TRYPSS	31A	06/11/2023
CONT1-OCFE07-TRYPSS	33B	17/01/2023
CONT 2-TRYPSS-US	33A	01/09/2022
BLISTER-3-WASH	26A	31/03/2025
TRYPSS-US MO	021	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B0/T %	72.19
BLANC		ng/mL	B/T %	1.27
CAL.1	36.00	ng/mL	B/B0	88.91
CAL.2	73.00	ng/mL	B/B0	78.97
CAL.3	160.00	ng/mL	B/B0	61.42
CAL.4	300.00	ng/mL	B/B0	46.24
CAL.5	630.00	ng/mL	B/B0	33.22
CAL.6	1 560.00	ng/mL	B/B0	25.24

Activité totale / Total activity : 37583 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	ng/mL	184.00 - 276.00	243.14
CONTROL 2	ng/mL	784.00 - 1176.00	862.14

Date du contrôle
Control Date

18/01/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

21/03/2022

