

# TRYPS-US

Protocole/Protocol : 3h + 30mn

Lot N°/Batch # : 374U

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCFE07-TRYPS	374A	11/07/2022
Ab-OCFE07-TRYPS	24A	07/12/2023
RIP-OCFE07-TRYPS	21A	05/03/2025
CAL0-OCFE07-TRYPS	31A	06/11/2023
CAL1-OCFE07-TRYPS	31A	06/11/2023
CAL2-OCFE07-TRYPS	31A	06/11/2023
CAL3-OCFE07-TRYPS	31A	06/11/2023
CAL4-OCFE07-TRYPS	31A	06/11/2023
CAL5-OCFE07-TRYPS	31A	06/11/2023
CAL6-OCFE07-TRYPS	31A	06/11/2023
CONT1-OCFE07-TRYPS	33B	17/01/2023
CONT 2-TRYPS-US	33B	10/02/2023
BLISTER-3-WASH	26A	31/03/2025
TRYPS-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 %	72.13
BLANC		ng/mL	B/T %	0.90
CAL.1	36.00	ng/mL	B/B0	89.88
CAL.2	73.00	ng/mL	B/B0	78.43
CAL.3	160.00	ng/mL	B/B0	61.20
CAL.4	300.00	ng/mL	B/B0	46.98
CAL.5	630.00	ng/mL	B/B0	34.53
CAL.6	1 560.00	ng/mL	B/B0	26.20

Activité totale / Total activity : 39434 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	ng/mL	184.00 - 276.00	241.76
CONTROL 2	ng/mL	784.00 - 1176.00	1 089.07

Date du contrôle  
Control Date

10/05/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**11/07/2022**

