

OCFE07-TRYPS

Protocole/Protocol : 3h + 30mn

Lot N°/Batch # : 374A

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
BLISTER-3-WASH	26A	31/03/2025
RG-TRYPSIN MO	016	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B0/T %	72.71
BLANC		ng/mL	B/T %	1.01
CAL.1	36.00	ng/mL	B/B0	89.41
CAL.2	73.00	ng/mL	B/B0	78.68
CAL.3	160.00	ng/mL	B/B0	61.15
CAL.4	300.00	ng/mL	B/B0	46.59
CAL.5	630.00	ng/mL	B/B0	33.94
CAL.6	1 560.00	ng/mL	B/B0	25.75

Activité totale / Total activity : 38115 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	184.00 - 276.00	232.70

Date du contrôle
Control Date

13/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

11/07/2022

