

GASK-PR

Protocole/Protocol : 2h

Lot N°/Batch # : 542A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-GASK-PR	542A	18/07/2022
Ab-GASK-PR	50A	09/12/2022
CAL0-GASK-PR	41A	19/03/2024
CAL1-GASK-PR	40A	21/09/2023
CAL2-GASK-PR	40A	21/09/2023
CAL3-GASK-PR	40A	21/09/2023
CAL4-GASK-PR	40A	21/09/2023
CAL5-GASK-PR	40A	21/09/2023
REAC-PR-100	81A	12/01/2023
GASK-PR MO	029	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	µU/mL	B0/T %	34.91
BLANC		µU/mL	B/T %	1.48
CAL.1	22.00	µU/mL	B/B0	93.46
CAL.2	60.00	µU/mL	B/B0	75.49
CAL.3	141.00	µU/mL	B/B0	55.14
CAL.4	460.00	µU/mL	B/B0	33.21
CAL.5	1 480.00	µU/mL	B/B0	16.47

Activité totale / Total activity : 18137 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
40391	µU/mL	36.00 - 70.00	62.67
40392	µU/mL	202.00 - 310.00	231.24
40393	µU/mL	410.00 - 630.00	546.46

Date du contrôle
Control Date

16/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

18/07/2022

