

GASK-PR

Protocole/Protocol : 2h

Lot N°/Batch # : 532A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-GASK-PR	532A	18/10/2021
Ab-GASK-PR	50A	09/12/2022
CAL0-GASK-PR	39A	04/07/2022
CAL1-GASK-PR	39B	19/01/2022
CAL2-GASK-PR	39B	19/01/2022
CAL3-GASK-PR	39B	19/01/2022
CAL4-GASK-PR	39B	19/01/2022
CAL5-GASK-PR	39B	19/01/2022
REAC-PR-100	80A	20/06/2022
GASK-PR MO	029	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
C.LIAISON	0.00	µU/mL	B0/T %	31.76
BLANC	0.00	µU/mL	B/T %	1.31
CAL.1	20.00	µU/mL	B/B0	96.87
CAL.2	54.00	µU/mL	B/B0	81.05
CAL.3	144.00	µU/mL	B/B0	58.38
CAL.4	480.00	µU/mL	B/B0	33.21
CAL.5	1 600.00	µU/mL	B/B0	17.86

Activité totale / Total activity : 18490 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
40391	µU/mL	36.00 - 70.00	60.07
40392	µU/mL	202.00 - 310.00	260.87
40393	µU/mL	410.00 - 630.00	578.52

Date du contrôle
Control Date

17/08/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

18/10/2021

