

GASK-PR

Protocole/Protocol : 2h

Lot N°/Batch # : 531A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-GASK-PR	531A	20/09/2021
Ab-GASK-PR	50A	09/12/2022
CAL0-GASK-PR	39A	04/07/2022
CAL1-GASK-PR	39B	19/01/2022
CAL2-GASK-PR	39C	19/01/2022
CAL3-GASK-PR	39C	19/01/2022
CAL4-GASK-PR	39C	19/01/2022
CAL5-GASK-PR	39C	19/01/2022
REAC-PR-100	80A	20/06/2022
GASK-PR MO	028	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
C.LIAISON	0.00	µU/mL	B0/T %	30.79
BLANC	0.00	µU/mL	B/T %	1.30
CAL.1	20.00	µU/mL	B/B0	98.03
CAL.2	50.00	µU/mL	B/B0	82.60
CAL.3	136.00	µU/mL	B/B0	59.79
CAL.4	450.00	µU/mL	B/B0	34.95
CAL.5	1 480.00	µU/mL	B/B0	17.62

Activité totale / Total activity : 18582 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
40391	µU/mL	36.00 - 70.00	56.10
40392	µU/mL	202.00 - 310.00	240.38
40393	µU/mL	410.00 - 630.00	523.24

Date du contrôle
Control Date

19/07/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

20/09/2021

