

RENINE

Protocole/Protocol : 3h

Lot N°/Batch # : 258A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-RENINE	258A	05/07/2022
CT-RENINE	82A	27/02/2024
CAL0-DIL-RENINE	19A	16/11/2022
CAL1-RENINE	25A	19/03/2024
CAL2-RENINE	25A	19/03/2024
CAL3-RENINE	25A	19/03/2024
CAL4-RENINE	25A	19/03/2024
CAL5-RENINE	25A	19/03/2024
CONT-RENINE	25A	19/03/2024
SOL-WASH-R3	16A	23/09/2022
RENIN III GENERATION	020	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	pg/mL	B/T %	0.02
CAL.1	2.70	pg/mL	B-B0/T	0.21
CAL.2	5.20	pg/mL	B-B0/T	0.45
CAL.3	21.70	pg/mL	B-B0/T	1.66
CAL.4	81.00	pg/mL	B-B0/T	6.59
CAL.5	280.00	pg/mL	B-B0/T	24.15

Activité totale / Total activity : 183302 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	18.90 - 28.30	20.95
AS008	pg/mL	3.00 - 4.60	3.52
DS007	pg/mL	62.20 - 93.20	76.22
ES007	pg/mL	114.40 - 171.60	141.21

Date du contrôle
Control Date

12/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

05/07/2022

