

RENINE

Protocole/Protocol : 3h

Lot N°/Batch # : 250A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-RENINE	250A	07/12/2021
CT-RENINE	81A	24/05/2022
CAL0-DIL-RENINE	19A	16/11/2022
CAL1-RENINE	22A	18/02/2022
CAL2-RENINE	22A	18/02/2022
CAL3-RENINE	22A	18/02/2022
CAL4-RENINE	22A	18/02/2022
CAL5-RENINE	22A	18/02/2022
CONT-RENINE	22A	18/02/2022
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 %	0.01
CAL.1	2.50	pg/mL	B-B0/T	0.17
CAL.2	5.00	pg/mL	B-B0/T	0.35
CAL.3	19.60	pg/mL	B-B0/T	1.41
CAL.4	78.00	pg/mL	B-B0/T	5.76
CAL.5	310.00	pg/mL	B-B0/T	24.53

Activité totale / Total activity : 183740 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	32.000 - 48.000	38.06
AS007	pg/mL	3.000 - 4.600	4.13
DS007	pg/mL	62.2 - 93.2	77.74

Date du contrôle
Control Date

14/10/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

07/12/2021

