

# RENINE

Protocole/Protocol : 3h

Lot N°/Batch # : 256B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-RENINE	256A	24/05/2022
CT-RENINE	82A	27/02/2024
CAL0-DIL-RENINE	19A	16/11/2022
CAL1-RENINE	23A	26/10/2023
CAL2-RENINE	23A	26/10/2023
CAL3-RENINE	23A	26/10/2023
CAL4-RENINE	23A	26/10/2023
CAL5-RENINE	23A	26/10/2023
CONT-RENINE	24A	01/12/2023
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	2.50	pg/mL	B-B0/T	0.20
CAL.2	5.00	pg/mL	B-B0/T	0.43
CAL.3	21.60	pg/mL	B-B0/T	1.80
CAL.4	88.00	pg/mL	B-B0/T	7.18
CAL.5	350.00	pg/mL	B-B0/T	29.49

Activité totale / Total activity : 176454 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	20.600 - 30.800	26.16
AS007	pg/mL	3.00 - 4.60	4.23
DS007	pg/mL	62.20 - 93.20	78.94
ES007	pg/mL	114.40 - 171.60	139.57

Date du contrôle  
Control Date

31/03/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

24/05/2022

