

RENINE

Protocole/Protocol : 3h

Lot N°/Batch # : 220A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-RENINE	220A	20/08/2019
CT-RENINE	78A	06/04/2021
CAL0-DIL-RENINE	18A	09/02/2021
CAL1-RENINE	21A	28/10/2020
CAL2-RENINE	21A	28/10/2020
CAL3-RENINE	21A	28/10/2020
CAL4-RENINE	21A	28/10/2020
CAL5-RENINE	21A	28/10/2020
CONT-RENINE	21A	28/10/2020
SOL-WASH-R3	15A	01/01/2021
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	2.80	pg/mL	B-B0/T	0.22
CAL.2	5.30	pg/mL	B-B0/T	0.45
CAL.3	20.60	pg/mL	B-B0/T	1.78
CAL.4	80.00	pg/mL	B-B0/T	7.07
CAL.5	310.00	pg/mL	B-B0/T	29.07

Activité totale / Total activity : 178312 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	29.600 - 44.400	36.81
AS005	pg/mL	3.00 - 4.60	4.28
ES006	pg/mL	124.80 - 187.20	158.67

Date du contrôle
Control Date

28/06/2019

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

20/08/2019

