

CORT-CT2

Protocole/Protocol : 2h

Lot N°/Batch # : 119A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-CORT-CT2	119A	04/08/2022
CT-CORT-CT2	20A	19/01/2023
CAL0-CORT-CT2	06A	12/01/2024
CAL1-CORT-CT2	06A	12/01/2024
CAL2-CORT-CT2	06A	12/01/2024
CAL3-CORT-CT2	06A	12/01/2024
CAL4-CORT-CT2	06A	12/01/2024
CORT CT2 RIA MO	025	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
Calibrator	0.00	nmol/L	B/T %	86.91
Calibrator 1	20.00	nmol/L	B/B0	84.33
Calibrator 2	73.00	nmol/L	B/B0	61.46
Calibrator 3	510.00	nmol/L	B/B0	30.35
Calibrator 4	2 000.00	nmol/L	B/B0	16.20

Activité totale / Total activity : 80460 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
OHC 1	nmol/L	52.10 - 71.80	68.44
OHC 2	nmol/L	390.00 - 544.00	502.03
OHC 3	nmol/L	636.00 - 961.00	820.24

Date du contrôle
Control Date

20/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

04/08/2022

