

TESTO-CT2

Protocole/Protocol : 1h

Lot N°/Batch # : 73B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TESTO CT2 TRACERS	73A	25/08/2022
CT-TESTO-CT2	34B	13/07/2023
CAL0-TESTO-CT2	08A	18/10/2023
CAL1-TESTO-CT2	08A	18/10/2023
CAL2-TESTO-CT2	08A	18/10/2023
CAL3-TESTO-CT2	08A	18/10/2023
CAL4-TESTO-CT2	08A	18/10/2023
CAL5-TESTO-CT2	08A	18/10/2023
TESTO CT2 RIA MO	026	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
Calibrator 0	0.00	nmol/L	B0/T %	43.54
Calibrator 1	0.40	nmol/L	B/B0	91.71
Calibrator 2	1.20	nmol/L	B/B0	80.12
Calibrator 3	4.80	nmol/L	B/B0	58.41
Calibrator 4	13.00	nmol/L	B/B0	39.54
Calibrator 5	43.00	nmol/L	B/B0	24.89

Activité totale / Total activity : 76915 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
OHC 1	nmol/L	0.64 - 1.28	1.09
OHC 2	nmol/L	4.07 - 5.85	4.74
OHC 3	nmol/L	29.40 - 40.20	34.30

Date du contrôle
Control Date

17/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

25/08/2022

