

# TESTO-CT2

Protocole/Protocol : 1h

Lot N°/Batch # : 73A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TESTO CT2 TRACERS	73A	25/08/2022
CT-TESTO-CT2	34A	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 %	44.47
Calibrator 1	0.40	nmol/L	B/B0	90.34
Calibrator 2	1.20	nmol/L	B/B0	79.41
Calibrator 3	4.80	nmol/L	B/B0	56.28
Calibrator 4	13.00	nmol/L	B/B0	36.93
Calibrator 5	43.00	nmol/L	B/B0	23.88

Activité totale / Total activity : 78052 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
OHC 1	nmol/L	0.64 - 1.28	1.10
OHC 2	nmol/L	4.07 - 5.85	5.10
OHC 3	nmol/L	29.40 - 40.20	33.50

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**

