

# TESTO-CT2

Protocole/Protocol : 1h

Lot N°/Batch # : 72A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TESTO CT2 TRACERS	72A	07/07/2022
CT-TESTO-CT2	33A	06/04/2023
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 %	38.13
Calibrator 1	0.40	nmol/L	B/B0	92.18
Calibrator 2	1.20	nmol/L	B/B0	80.98
Calibrator 3	4.80	nmol/L	B/B0	61.05
Calibrator 4	13.00	nmol/L	B/B0	43.49
Calibrator 5	43.00	nmol/L	B/B0	26.80

Activité totale / Total activity : 84222 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
OHC 1 F356A	nmol/L	0.64 - 1.28	0.88
OHC 2 F356B	nmol/L	4.07 - 5.85	4.52
OHC 3 F356C	nmol/L	29.40 - 40.20	32.52

Date du contrôle  
Control Date

28/03/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

07/07/2022

