

# TESTO-CT2

Protocole/Protocol : 1h

Lot N°/Batch # : 69A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TESTO CT2 TRACERS	69A	20/01/2022
CT-TESTO-CT2	32A	05/01/2023
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 %	40.32
Calibrator 1	0.40	nmol/L	B/B0	91.61
Calibrator 2	1.20	nmol/L	B/B0	81.74
Calibrator 3	4.80	nmol/L	B/B0	59.76
Calibrator 4	13.00	nmol/L	B/B0	39.30
Calibrator 5	43.00	nmol/L	B/B0	26.14

Activité totale / Total activity : 76559 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
OHC 1	nmol/L	0.64 - 1.28	1.03
OHC 2	nmol/L	4.07 - 5.85	5.32
OHC 3	nmol/L	29.40 - 40.20	31.00

Date du contrôle  
Control Date

12/10/2021

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

20/01/2022

