

TESTO-CT2

Protocole/Protocol : 1h

Lot N°/Batch # : 67A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TESTO CT2 TRACERS	67A	30/09/2021
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 %	38.65
Calibrator 1	0.40	nmol/L	B/B0	91.08
Calibrator 2	1.20	nmol/L	B/B0	81.58
Calibrator 3	4.80	nmol/L	B/B0	60.32
Calibrator 4	13.00	nmol/L	B/B0	40.61
Calibrator 5	43.00	nmol/L	B/B0	26.59

Activité totale / Total activity : 78853 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
OHC 1	nmol/L	0.64 - 1.28	0.85
OHC 2	nmol/L	4.07 - 5.85	4.84
OHC 3	nmol/L	29.40 - 40.20	31.89

Date du contrôle
Control Date

18/06/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

30/09/2021

