

ESTR-US-CT

Protocole/Protocol : 1h + 1h

Lot N°/Batch # : 113A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
ESTR US CT TRACER	113A	09/12/2021
CT-ESTR-US-CT	38A	26/04/2022
BUF-ESTR-US-CT	19A	03/03/2022
CAL0-ESTR-US-CT	17A	30/09/2022
CAL1-ESTR-US-CT	17A	30/09/2022
CAL2-ESTR-US-CT	18B	04/03/2023
CAL3-ESTR-US-CT	17A	30/09/2022
CAL4-ESTR-US-CT	17A	30/09/2022
CAL5-ESTR-US-CT	17A	30/09/2022
TWEEN-1-3-EL-1FLA	215A	03/06/2024
ESTR-US-CT MO	021	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
Calibrator	0.00	pmol/l	B/T %	36.44
Calibrator 1	12.00	pmol/l	B/B0	92.47
Calibrator 2	30.00	pmol/l	B/B0	79.71
Calibrator 3	90.00	pmol/l	B/B0	60.14
Calibrator 4	400.00	pmol/l	B/B0	29.85
Calibrator 5	1 890.00	pmol/l	B/B0	10.55

Activité totale / Total activity : 17115 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
40371	pmol/L	138.00 - 208.00	187.14
40372	pmol/L	310.00 - 470.00	379.56
40373	pmol/L	480.00 - 720.00	692.20

Date du contrôle
Control Date

08/10/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

09/12/2021

