

THYRO

Protocole/Protocol : 3h + 16 - 20h

Lot N°/Batch # : 255A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-THYRO	255A	07/03/2022
CT-THYRO	80A	06/03/2023
BUF-THYRO	25A	27/06/2022
CAL1-THYRO	36G	24/03/2022
CAL2-THYRO	36G	24/03/2022
CAL3-THYRO	36G	24/03/2022
CAL4-THYRO	36G	24/03/2022
CAL5-THYRO	36G	24/03/2022
CAL6-THYRO	36G	24/03/2022
CAL7-THYRO	36G	24/03/2022
CONT1-THYRO	18F	10/11/2022
CONT2-THYRO	18F	10/11/2022
SOL-WASH-R3	16A	23/09/2022
THYRO IRMA MO	018	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.1	0.21	ng/mL	B/T %	0.14
CAL.2	1.95	ng/mL	B/T %	0.34
CAL.3	5.20	ng/mL	B/T %	0.75
CAL.4	16.00	ng/mL	B/T %	2.00
CAL.5	56.00	ng/mL	B/T %	6.84
CAL.6	200.00	ng/mL	B/T %	26.87
CAL.7	480.00	ng/mL	B/T %	55.35

Activité totale / Total activity : 254423 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	ng/mL	9.92 - 14.88	13.59
CONTROL 2	ng/mL	32.00 - 48.00	42.89

Date du contrôle
Control Date

10/01/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

07/03/2022

