

THYRO

Protocole/Protocol : 3h + 16 - 20h

Lot N°/Batch # : 249B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-THYRO	249A	20/09/2021
CT-THYRO	79A	12/12/2022
BUF-THYRO	25A	27/06/2022
CAL1-THYRO	36B	24/03/2022
CAL2-THYRO	36B	24/03/2022
CAL3-THYRO	36B	24/03/2022
CAL4-THYRO	36B	24/03/2022
CAL5-THYRO	36B	24/03/2022
CAL6-THYRO	36B	24/03/2022
CAL7-THYRO	36B	24/03/2022
CONT1-THYRO	18B	10/11/2022
CONT2-THYRO	18C	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 %	0.13
CAL.2	1.75	ng/mL	B/T %	0.31
CAL.3	5.50	ng/mL	B/T %	0.70
CAL.4	17.00	ng/mL	B/T %	1.92
CAL.5	61.00	ng/mL	B/T %	6.66
CAL.6	220.00	ng/mL	B/T %	25.91
CAL.7	480.00	ng/mL	B/T %	53.94

Activité totale / Total activity : 254761 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	ng/mL	11.52 - 17.28	16.00
CONTROL 2	ng/mL	35.20 - 52.80	50.34

Date du contrôle
Control Date

27/07/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

20/09/2021

