

THYRO

Protocole/Protocol : 3h + 16 - 20h

Lot N°/Batch # : 259A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-THYRO	259A	04/07/2022
CT-THYRO	81A	06/06/2023
BUF-THYRO	26A	26/09/2023
CAL1-THYRO	37A	02/12/2022
CAL2-THYRO	37A	02/12/2022
CAL3-THYRO	37B	02/12/2022
CAL4-THYRO	37B	02/12/2022
CAL5-THYRO	37A	02/12/2022
CAL6-THYRO	37A	02/12/2022
CAL7-THYRO	37A	02/12/2022
CONT1-THYRO	18G	10/11/2022
CONT2-THYRO	18G	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 %	0.17
CAL.2	1.67	ng/mL	B/T %	0.32
CAL.3	5.50	ng/mL	B/T %	0.77
CAL.4	14.60	ng/mL	B/T %	1.66
CAL.5	51.00	ng/mL	B/T %	6.11
CAL.6	196.00	ng/mL	B/T %	24.45
CAL.7	520.00	ng/mL	B/T %	55.35

Activité totale / Total activity : 243818 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	ng/mL	9.52 - 14.28	13.57
CONTROL 2	ng/mL	29.60 - 44.40	39.66

Date du contrôle
Control Date

06/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

04/07/2022

