

ELSA-FBHCG

Protocole/Protocol : 1h + 2h

Lot N°/Batch # : 435A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-FBHCG	435A	07/07/2022
CT-ELSA-FBHCG	110F	08/08/2022
BUF-ELSA-FBHCG	57A	05/01/2023
CAL0-ELSA-FBHCG	69A	10/01/2023
CAL1-ELSA-FBHCG	68A	12/01/2023
CAL2-ELSA-FBHCG	68A	12/01/2023
CAL3-ELSA-FBHCG	68A	12/01/2023
CAL4-ELSA-FBHCG	68A	12/01/2023
CAL5-ELSA-FBHCG	68A	12/01/2023
CONT-ELSA-FBHCG	68A	12/01/2023
TWEEN-1-3-EL-1FLA	215A	03/06/2024
ELSA-F-BHCG MO	023	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B/T %	0.06
CAL.1	0.22	ng/mL	B/T %	0.42
CAL.2	2.10	ng/mL	B/T %	4.08
CAL.3	9.00	ng/mL	B/T %	17.16
CAL.4	22.00	ng/mL	B/T %	36.91
CAL.5	50.00	ng/mL	B/T %	63.38

Activité totale / Total activity : 274389 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	0.91 - 1.37	1.10

Date du contrôle
Control Date

10/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

07/07/2022

