

CGA-ELISA

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 75A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA	75A	19/09/2023
MICROPLATE-CGA-ELISA	80A	20/09/2023
CAL0-DIL-CGA-ELISA	39A	12/09/2023
BUF-CONJ-CGA-ELISA	26A	10/08/2023
CAL1-CGA-ELISA	23A	23/01/2024
CAL2-CGA-ELISA	23A	23/01/2024
CAL3-CGA-ELISA	23A	23/01/2024
CAL4-CGA-ELISA	23A	23/01/2024
CAL5-CGA-ELISA	23A	23/01/2024
CONT-CGA-ELISA	23A	23/01/2024
CONT2-CGA-ELISA	23A	23/01/2024
TWEEN-1-3-EL-1FLA	215A	03/06/2024
SUBS-TMB	57A	29/10/2024
STOP-SOL	31A	28/12/2023
CHROMO@(CGA-ELISA)MO	017	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL0	0.00	ng/mL	DO	0.04
CAL1	61.00	ng/mL	DO	0.16
CAL2	135.00	ng/mL	DO	0.38
CAL3	270.00	ng/mL	DO	0.74
CAL4	490.00	ng/mL	DO	1.46
CAL5	850.00	ng/mL	DO	2.36

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	72 - 108	92.87
CONTROL2	ng/mL	496 - 744	604.80

Date du contrôle
Control Date

16/02/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

10/08/2023

