

CGA-ELISA

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 77A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA	77A	12/02/2024
MICROPLATE-CGA-ELISA	81A	01/02/2024
CAL0-DIL-CGA-ELISA	40A	13/02/2024
BUF-CONJ-CGA-ELISA	27A	27/07/2024
CAL1-CGA-ELISA	24A	14/06/2024
CAL2-CGA-ELISA	24A	14/06/2024
CAL3-CGA-ELISA	24A	14/06/2024
CAL4-CGA-ELISA	24A	14/06/2024
CAL5-CGA-ELISA	24A	14/06/2024
CONT-CGA-ELISA	24A	14/06/2024
CONT2-CGA-ELISA	24A	14/06/2024
TWEEN 20	215A	03/06/2024
SUBS TMB	59A	27/01/2025
STOP SOLN	33A	05/06/2024
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	0.03
CAL1	76.00	ng/mL	DO	0.18
CAL2	145.00	ng/mL	DO	0.36
CAL3	240.00	ng/mL	DO	0.63
CAL4	480.00	ng/mL	DO	1.31
CAL5	850.00	ng/mL	DO	2.16

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	81.6 - 122.4	100.34
CONTROL2	ng/mL	480 - 720	594.66

Date du contrôle
Control Date

18/07/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

01/02/2024

