

CGA-ELISA

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 72A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA	72A	14/02/2023
MICROPLATE-CGA-ELISA	79A	09/02/2023
CAL0-DIL-CGA-ELISA	38A	13/02/2023
BUF-CONJ-CGA-ELISA	25A	10/08/2023
CAL1-CGA-ELISA	22A	29/05/2023
CAL2-CGA-ELISA	22A	29/05/2023
CAL3-CGA-ELISA	22A	29/05/2023
CAL4-CGA-ELISA	22A	29/05/2023
CAL5-CGA-ELISA	22A	29/05/2023
CONT-CGA-ELISA	22A	29/05/2023
CONT2-CGA-ELISA	22A	29/05/2023
TWEEN-1-3-EL-1FLA	215A	03/06/2024
SUBS-TMB	56A	17/12/2023
STOP-SOL	30A	18/05/2023
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	0.03
CAL1	62.00	ng/mL	DO	0.19
CAL2	120.00	ng/mL	DO	0.40
CAL3	220.00	ng/mL	DO	0.79
CAL4	430.00	ng/mL	DO	1.59
CAL5	760.00	ng/mL	DO	2.67

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	68.8 - 103.2	81.81
CONTROL2	ng/mL	424 - 636	522.98

Date du contrôle
Control Date

13/07/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

09/02/2023

