

# CGA-ELISA

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 71A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA	71A	21/11/2022
MICROPLATE-CGA-ELISA	78A	07/07/2022
CAL0-DIL-CGA-ELISA	37A	04/07/2022
BUF-CONJ-CGA-ELISA	24A	06/08/2022
CAL1-CGA-ELISA	21B	06/10/2022
CAL2-CGA-ELISA	21B	06/10/2022
CAL3-CGA-ELISA	21B	06/10/2022
CAL4-CGA-ELISA	21B	06/10/2022
CAL5-CGA-ELISA	21B	06/10/2022
CONT-CGA-ELISA	21B	06/10/2022
CONT2-CGA-ELISA	21B	06/10/2022
TWEEN-1-3-EL-1FLA	213A	20/03/2023
SUBS-TMB	55A	08/07/2023
STOP-SOL	29A	01/11/2022
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	60.00	ng/mL	DO	0.14
CAL2	130.00	ng/mL	DO	0.33
CAL3	240.00	ng/mL	DO	0.63
CAL4	460.00	ng/mL	DO	1.30
CAL5	850.00	ng/mL	DO	2.41

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	65.6 - 98.4	89.13
CONTROL2	ng/mL	480 - 720	585.33

Date du contrôle  
Control Date

05/05/2021

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**04/07/2022**

