

CGA-ELISA

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 70A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA	70A	09/08/2022
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	21A	06/10/2022
CAL2-CGA-ELISA	21A	06/10/2022
CAL3-CGA-ELISA	21A	06/10/2022
CAL4-CGA-ELISA	21A	06/10/2022
CAL5-CGA-ELISA	21A	06/10/2022
CONT-CGA-ELISA	21A	06/10/2022
CONT2-CGA-ELISA	21A	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	0.02
CAL1	60.00	ng/mL	DO	0.14
CAL2	130.00	ng/mL	DO	0.36
CAL3	220.00	ng/mL	DO	0.67
CAL4	480.00	ng/mL	DO	1.66
CAL5	900.00	ng/mL	DO	2.55

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	70.400 - 105.600	92.64
CONTROL2	ng/mL	432 - 648	512.82

Date du contrôle
Control Date

18/12/2020

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

04/07/2022

