

# CGA-ELISA

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 60C

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA	60A	07/10/2021
MICROPLATE-CGA-ELISA	71A	17/11/2021
CAL0-DIL-CGA-ELISA	36A	23/12/2021
BUF-CONJ-CGA-ELISA	23A	06/12/2021
CAL1-CGA-ELISA	18A	01/10/2021
CAL2-CGA-ELISA	18A	01/10/2021
CAL3-CGA-ELISA	18A	01/10/2021
CAL4-CGA-ELISA	18A	01/10/2021
CAL5-CGA-ELISA	18A	01/10/2021
CONT-CGA-ELISA	18A	01/10/2021
CONT2-CGA-ELISA	18A	01/10/2021
TWEEN-1-3-EL-1FLA	213A	20/03/2023
STOP-SOL	28A	28/03/2022
SUBS-TMB	54A	27/02/2023
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	0.01
CAL1	70.00	ng/mL	DO	0.13
CAL2	135.00	ng/mL	DO	0.29
CAL3	230.00	ng/mL	DO	0.60
CAL4	410.00	ng/mL	DO	1.16
CAL5	720.00	ng/mL	DO	2.14

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	80 - 120	93.60
CONTROL2	ng/mL	392 - 588	504.33

Date du contrôle  
Control Date

17/06/2020

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

01/10/2021

