

CGA-ELISA

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 66A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA	66A	16/05/2022
MICROPLATE-CGA-ELISA	75A	31/03/2022
CAL0-DIL-CGA-ELISA	36A	23/12/2021
BUF-CONJ-CGA-ELISA	24A	06/08/2022
CAL1-CGA-ELISA	19A	18/04/2022
CAL2-CGA-ELISA	19A	18/04/2022
CAL3-CGA-ELISA	19A	18/04/2022
CAL4-CGA-ELISA	19A	18/04/2022
CAL5-CGA-ELISA	19A	18/04/2022
CONT-CGA-ELISA	19A	18/04/2022
CONT2-CGA-ELISA	19A	18/04/2022
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	62.00	ng/mL	DO	0.16
CAL2	130.00	ng/mL	DO	0.34
CAL3	240.00	ng/mL	DO	0.65
CAL4	500.00	ng/mL	DO	1.37
CAL5	900.00	ng/mL	DO	2.53

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	65.6 - 98.4	76.92
CONTROL2	ng/mL	528 - 792	674.56

Date du contrôle
Control Date

29/09/2020

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

23/12/2021

