

CGA-ELISA

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 60B

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	35A	16/07/2021
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	211A	05/01/2023
STOP-SOL	27A	14/10/2021
SUBS-TMB	52A	22/05/2022
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	0.01
CAL1	70.00	ng/mL	DO	0.15
CAL2	135.00	ng/mL	DO	0.33
CAL3	230.00	ng/mL	DO	0.66
CAL4	410.00	ng/mL	DO	1.29
CAL5	720.00	ng/mL	DO	2.05

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	80 - 120	106.28
CONTROL2	ng/mL	392 - 588	498.79

Date du contrôle
Control Date

14/04/2020

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

16/07/2021

