

CGA-ELISA-US

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 68W

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA-US	68V	12/07/2022
MICROPLAQUE-CGA-ELIS	77V	30/06/2022
CAL0-DIL-CGA-ELISA-U	37V	04/07/2022
CAL1-CGA-ELISA-US	21V	06/10/2022
CAL2-CGA-ELISA-US	21V	06/10/2022
CAL3-CGA-ELISA-US	21V	06/10/2022
CAL4-CGA-ELISA-US	21V	06/10/2022
CAL5-CGA-ELISA-US	21V	06/10/2022
CONT1-CGA-ELISA-US	21V	06/10/2022
CONT2-CGA-ELISA-US	21V	06/10/2022
TWEEN-20-US	213V	20/03/2023
SUBS-TMB-US	55V	08/07/2023
STOP-SOL-US	29V	01/11/2022
CGA elisa MO	017	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL0	0.00	ng/mL	DO	0.05
CAL1	47.00	ng/mL	DO	0.13
CAL2	100.00	ng/mL	DO	0.28
CAL3	193.00	ng/mL	DO	0.65
CAL4	360.00	ng/mL	DO	1.17
CAL5	770.00	ng/mL	DO	2.27

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	56 - 84	70.15
CONTROL 2	ng/mL	376 - 564	490.18

Date du contrôle
Control Date

14/12/2020

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

30/06/2022

