

CGA-ELISA-US

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 73V

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA-US	73V	13/02/2023
MICROPLAQUE-CGA-ELIS	79V	09/02/2023
CAL0-DIL-CGA-ELISA-U	38V	13/02/2023
CAL1-CGA-ELISA-US	22V	29/05/2023
CAL2-CGA-ELISA-US	22V	29/05/2023
CAL3-CGA-ELISA-US	22V	29/05/2023
CAL4-CGA-ELISA-US	22V	29/05/2023
CAL5-CGA-ELISA-US	22V	29/05/2023
CONT1-CGA-ELISA-US	22V	29/05/2023
CONT2-CGA-ELISA-US	22V	29/05/2023
TWEEN-20-US	213V	20/03/2023
SUBS-TMB-US	55V	08/07/2023
STOP-SOL-US	30V	18/05/2023
CGA elisa MO	018	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL0	0.00	ng/mL	DO	0.05
CAL1	61.00	ng/mL	DO	0.18
CAL2	130.00	ng/mL	DO	0.36
CAL3	230.00	ng/mL	DO	0.67
CAL4	450.00	ng/mL	DO	1.36
CAL5	760.00	ng/mL	DO	2.41

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	68.8 - 103.2	84.77
CONTROL 2	ng/mL	440 - 660	551.64

Date du contrôle
Control Date

20/07/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

09/02/2023

