

CGA-ELISA-US

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 78V

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA-US	78V	06/02/2024
MICROPLAQUE-CGA-ELIS	81V	01/02/2024
CAL0-DIL-CGA-ELISA-U	40V	13/02/2024
CAL1-CGA-ELISA-US	24V	14/06/2024
CAL2-CGA-ELISA-US	24V	14/06/2024
CAL3-CGA-ELISA-US	24V	14/06/2024
CAL4-CGA-ELISA-US	24V	14/06/2024
CAL5-CGA-ELISA-US	24V	14/06/2024
CONT1-CGA-ELISA-US	24V	14/06/2024
CONT2-CGA-ELISA-US	24V	14/06/2024
TWEEN 20	215V	03/06/2024
SUBS TMB	58V	29/10/2024
STOP SOLN	29V	30/04/2024
CGA elisa MO	018	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL0	0.00	ng/mL	DO	0.07
CAL1	59.00	ng/mL	DO	0.18
CAL2	118.00	ng/mL	DO	0.34
CAL3	207.00	ng/mL	DO	0.58
CAL4	400.00	ng/mL	DO	1.20
CAL5	740.00	ng/mL	DO	2.15

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	64 - 96	77.82
CONTROL 2	ng/mL	416 - 624	516.00

Date du contrôle
Control Date

18/07/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

01/02/2024



DELEUZE Isabelle