

CGA-ELISA-NG

Protocole/Protocol : 1H + 2H

Lot N°/Batch # : 04B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA-NG	04A	11/04/2024
MICROPLATE-CGA-ELISA	04A	04/04/2024
DIL CAL0-CGA-ELISA-N	41NG	05/04/2024
CAL1-CGA-ELISA-NG	04A	11/04/2024
CAL2-CGA-ELISA-NG	04A	11/04/2024
CAL3-CGA-ELISA-NG	04A	11/04/2024
CAL4-CGA-ELISA-NG	04A	11/04/2024
CAL5-CGA-ELISA-NG	04A	11/04/2024
CONT1-CGA-ELISA-NG	04A	11/04/2024
CONT2-CGA-ELISA-NG	04A	11/04/2024
TWEEN-1-3-EL-1FLA	215AP	03/06/2024
SUBS-TMB	60AP	24/06/2024
STOP-SOLN	34AP	08/08/2024
CGA-ELISA-NG MO	03	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL0	0.00	ng/mL	DO	0.10
CAL1	80.00	ng/mL	DO	0.18
CAL2	160.00	ng/mL	DO	0.31
CAL3	320.00	ng/mL	DO	0.58
CAL4	600.00	ng/mL	DO	1.13
CAL5	1 100.00	ng/mL	DO	1.93

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	80.000 - 120.000	97.36
CONTROL2	ng/mL	608.000 - 912.000	753.70

Date du contrôle
Control Date

02/01/2023

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

04/04/2024

