

CGA-ELISA-NG

Protocole/Protocol : 1H + 2H

Lot N°/Batch # : 05B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ-CGA-ELISA-NG	05A	04/10/2024
MICROPLATE-CGA-ELISA	05A	02/10/2024
DIL CAL0-CGA-ELISA-N	42A	10/10/2024
CAL1-CGA-ELISA-NG	05A	10/10/2024
CAL2-CGA-ELISA-NG	05A	10/10/2024
CAL3-CGA-ELISA-NG	05A	10/10/2024
CAL4-CGA-ELISA-NG	05A	10/10/2024
CAL5-CGA-ELISA-NG	05A	10/10/2024
CONT1-CGA-ELISA-NG	05A	10/10/2024
CONT2-CGA-ELISA-NG	05A	10/10/2024
TWEEN 20	216A	18/06/2026
STOP SOLN	35AP	07/02/2025
SUBS TMB	61AP	27/10/2025
CGA-ELISA-NG MO	004	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL0	0.00	ng/mL	DO	0.09
CAL1	54.00	ng/mL	DO	0.16
CAL2	118.00	ng/mL	DO	0.28
CAL3	293.00	ng/mL	DO	0.66
CAL4	610.00	ng/mL	DO	1.36
CAL5	1 010.00	ng/mL	DO	2.28

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	59.200 - 88.800	73.16
CONTROL2	ng/mL	592.000 - 888.000	729.46

Date du contrôle
Control Date

06/09/2023

Date de péremption
Expiry date

02/10/2024

Visa du responsable du contrôle qualité
Quality control manager visa



Cisbio Bioassays
Parc Marcel Boiteux - BP 84175
30200 Codolet / France



DELEUZE MAZZOCCA ISABELLE