

In-vitro Diagnostic Reagents

CGA-ELISA-NG

Product number:	CGA-ELISA-NG	Lot number:	06B	Protocol:	1H + 2H
Reagents	Batch #	Expiry Date			
Réactifs	Lot N°	Péremption			
CONJ-CGA-ELISA-NG	06B	18/08/2025			
MICROPLATE-CGA-ELISA	06A	19/08/2025			
DIL CAL0-CGA-ELISA-N	43B	31/07/2025			
CAL1-CGA-ELISA-NG	06B	04/08/2025			
CAL2-CGA-ELISA-NG	06B	04/08/2025			
CAL3-CGA-ELISA-NG	06B	04/08/2025			
CAL4-CGA-ELISA-NG	06B	04/08/2025			
CAL5-CGA-ELISA-NG	06B	04/08/2025			
CONT1-CGA-ELISA-NG	06B	04/08/2025			
CONT2-CGA-ELISA-NG	06B	04/08/2025			
TWEEN 20	216B	18/06/2026			
SUBS TMB	62B	08/12/2025			
STOP SOLN	36B	01/01/2026			
CGA-ELISA-NG MO	05				

Results / Résultats

Parameter	Conc.	Unit	Response	Measured value
Paramètre	Conc.	Unité	Réponse	Valeur mesurée
CAL0	0.00	ng/mL	DO	0.11
CAL1	95.00	ng/mL	DO	0.23
CAL2	185.00	ng/mL	DO	0.41
CAL3	370.00	ng/mL	DO	0.83
CAL4	740.00	ng/mL	DO	1.68
CAL5	1 330.00	ng/mL	DO	2.57

Controls	Unit	Expected values	Measured value
Contrôles	Unité	Valeurs attendues	Valeur mesurée
CONTROL	ng/mL	96.000 - 144.000	118.82
CONTROL2	ng/mL	696.000 - 1044.000	848.88

Control Date 16/01/2024

Date du contrôle

Expiry date 31/07/2025

Date de péremption

Quality control manager visa

Visa du responsable du contrôle qualité



DELEUZE MAZZOCCA Isabelle