

In-vitro Diagnostic Reagents

CGA-ELISA-NG

Product number:	CGA-ELISA-NG	Lot number:	06A	Protocol:	1H + 2H
Reagents	Batch #			Expiry Date	
Réactifs	Lot N°			Péremption	
CONJ-CGA-ELISA-NG	06A			18/08/2025	
MICROPLATE-CGA-ELISA	06A			19/08/2025	
DIL CAL0-CGA-ELISA-N	43A			31/07/2025	
CAL1-CGA-ELISA-NG	06A			04/08/2025	
CAL2-CGA-ELISA-NG	06A			04/08/2025	
CAL3-CGA-ELISA-NG	06A			04/08/2025	
CAL4-CGA-ELISA-NG	06A			04/08/2025	
CAL5-CGA-ELISA-NG	06A			04/08/2025	
CONT1-CGA-ELISA-NG	06A			04/08/2025	
CONT2-CGA-ELISA-NG	06A			04/08/2025	
TWEEN 20	216A			18/06/2026	
SUBS TMB	62A			08/12/2025	
STOP SOLN	36A			01/01/2026	
CGA-ELISA-NG MO	004				

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	0.22
CAL2	185.00	ng/mL	DO	0.38
CAL3	370.00	ng/mL	DO	0.80
CAL4	740.00	ng/mL	DO	1.44
CAL5	1 330.00	ng/mL	DO	2.41

Controls	Unit	Expected values	Measured value
Contrôles	Unité	Valeurs attendues	Valeur mesurée
CONTROL	ng/mL	96.000 - 144.000	123.70
CONTROL2	ng/mL	696.000 - 1044.000	853.43

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