

In-vitro Diagnostic Reagents

CGA-ELISA-NG

Product number:	CGA-ELISA-NG	Lot number:	05C	Protocol:	1H + 2H
Reagents	Batch #	Expiry Date			
Réactifs	Lot N°	Péremption			
CONJ-CGA-ELISA-NG	05B	04/10/2024			
MICROPLATE-CGA-ELISA	05A	02/10/2024			
DIL CAL0-CGA-ELISA-N	42B	10/10/2024			
CAL1-CGA-ELISA-NG	05B	10/10/2024			
CAL2-CGA-ELISA-NG	05B	10/10/2024			
CAL3-CGA-ELISA-NG	05B	10/10/2024			
CAL4-CGA-ELISA-NG	05B	10/10/2024			
CAL5-CGA-ELISA-NG	05B	10/10/2024			
CONT1-CGA-ELISA-NG	05B	10/10/2024			
CONT2-CGA-ELISA-NG	05B	10/10/2024			
TWEEN 20	216A	18/06/2026			
SUBS TMB	61AP	27/10/2025			
STOP SOLN	35AP	07/02/2025			
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	0.09
CAL1	54.00	ng/mL	DO	0.17
CAL2	118.00	ng/mL	DO	0.29
CAL3	293.00	ng/mL	DO	0.72
CAL4	610.00	ng/mL	DO	1.43
CAL5	1 010.00	ng/mL	DO	2.43

Controls	Unit	Expected values	Measured value
Contrôles	Unité	Valeurs attendues	Valeur mesurée
CONTROL	ng/mL	59.2 - 88.8	80.22
CONTROL2	ng/mL	592 - 888	751.17

Control Date 24/01/2024

Date du contrôle

Quality control manager visa

Visa du responsable du contrôle qualité

Expiry date 02/10/2024

Date de péremption



DELEUZE MAZZOCCA Isabelle