

# CGA-ELISA-NG-US

Protocole/Protocol : 1H + 2H

Lot N°/Batch # : 04X

| Réactifs<br>Reagents | Lot N°<br>Batch # | Péremption<br>Expiry Date |
|----------------------|-------------------|---------------------------|
| CONJ-CGA-ELISA-NG-US | 04X               | 11/04/2024                |
| MICROPLATE-CGA-ELISA | 04X               | 04/04/2024                |
| DIL CAL0-CGA-ELISA-N | 41NGX             | 05/04/2024                |
| CAL1-CGA-ELISA-NG-US | 04X               | 11/04/2024                |
| CAL2-CGA-ELISA-NG-US | 04X               | 11/04/2024                |
| CAL3-CGA-ELISA-NG-US | 04X               | 11/04/2024                |
| CAL4-CGA-ELISA-NG-US | 04X               | 11/04/2024                |
| CAL5-CGA-ELISA-NG-US | 04X               | 11/04/2024                |
| CONT1-CGA-ELISA-NG-U | 04X               | 11/04/2024                |
| CONT2-CGA-ELISA-NG-U | 04X               | 11/04/2024                |
| TWEEN 20             | 215VP             | 03/06/2024                |
| SUBS TMB             | 60VP              | 24/06/2024                |
| STOP SOLN            | 34VP              | 08/08/2024                |
| CGA-ELISA-NG-US MO   | 01                |                           |

## RESULTATS / RESULTS

| Paramètre<br>Parameter | Conc.<br>Conc. | Unité<br>Unit | Réponse<br>Response | Valeur mesurée<br>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<br>Controls | Unité<br>Unit | Valeurs attendues<br>Expected values | Valeur mesurée<br>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

21/09/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

04/04/2024

