

# CGA-RIACT

Protocole/Protocol : 18 - 20h + 2h

Lot N°/Batch # : 211B

| Réactifs<br>Reagents | Lot N°<br>Batch # | Péremption<br>Expiry Date |
|----------------------|-------------------|---------------------------|
| TRA-CGA-RIACT        | 211A              | 01/07/2022                |
| CT-CGA-RIACT         | 48A               | 08/03/2023                |
| BUF-CGA-RIACT        | 45V               | 01/07/2022                |
| CAL1-CGA-RIACT       | 26A               | 16/02/2023                |
| CAL2-CGA-RIACT       | 26A               | 16/02/2023                |
| CAL3-CGA-RIACT       | 26A               | 16/02/2023                |
| CAL4-CGA-RIACT       | 26A               | 16/02/2023                |
| CAL5-CGA-RIACT       | 26A               | 16/02/2023                |
| CONT-CGA-RIACT       | 26A               | 16/02/2023                |
| BLISTER-3-WASH       | 26A               | 31/03/2025                |
| CGA-RIA-CT MO        | 026               |                           |

## RESULTATS / RESULTS

| Paramètre<br>Parameter | Conc.<br>Conc. | Unité<br>Unit | Réponse<br>Response | Valeur mesurée<br>Measured value |
|------------------------|----------------|---------------|---------------------|----------------------------------|
| CAL.0                  | 0.00           | ng/mL         | B/T %               | 0.02                             |
| CAL.1                  | 43.00          | ng/mL         | B/T %               | 2.21                             |
| CAL.2                  | 105.00         | ng/mL         | B/T %               | 5.85                             |
| CAL.3                  | 264.00         | ng/mL         | B/T %               | 15.69                            |
| CAL.4                  | 560.00         | ng/mL         | B/T %               | 31.47                            |
| CAL.5                  | 1 050.00       | ng/mL         | B/T %               | 59.58                            |

Activité totale / Total activity : 187415 cpm

| Contrôles<br>Controls | Unité<br>Unit | Valeurs attendues<br>Expected values | Valeur mesurée<br>Measured value |
|-----------------------|---------------|--------------------------------------|----------------------------------|
| CONTROL               | ng/mL         | 124.000 - 186.000                    | 156.71                           |

Date du contrôle  
Control Date

13/04/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

01/07/2022

