

CGA-RIACT

Protocole/Protocol : 18 - 20h + 2h

Lot N°/Batch # : 208A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-CGA-RIACT	208A	14/01/2022
CT-CGA-RIACT	49A	20/09/2023
BUF-CGA-RIACT	45A	13/06/2022
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-5-WASH	25B	28/02/2022
CGA-RIA-CT MO	025	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B/T %	0.02
CAL.1	43.00	ng/mL	B/T %	2.24
CAL.2	105.00	ng/mL	B/T %	5.79
CAL.3	264.00	ng/mL	B/T %	14.71
CAL.4	560.00	ng/mL	B/T %	30.68
CAL.5	1 050.00	ng/mL	B/T %	59.62

Activité totale / Total activity : 172401 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	124.000 - 186.000	155.62

Date du contrôle
Control Date

03/11/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

14/01/2022



DELEUZE Isabelle