

CEA-RIACT

Protocole/Protocol : 1h

Lot N°/Batch # : 497A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-CEA-RIACT	497A	04/07/2022
CT-CEA-RIACT	62A	18/01/2023
DIL-CEA-RIACT	28A	24/08/2022
CAL0-CEA-RIACT	30A	13/01/2023
CAL1-CEA-RIACT	30A	13/01/2023
CAL2-CEA-RIACT	30A	13/01/2023
CAL3-CEA-RIACT	30A	13/01/2023
CAL4-CEA-RIACT	30A	13/01/2023
CAL5-CEA-RIACT	30A	13/01/2023
CONT-CEA-RIACT	30A	13/01/2023
TWEEN-1-3-EL-1FLA	215A	03/06/2024
CEA-RIACT MO	012	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B/T %	0.12
CAL.1	2.70	ng/mL	B/T %	0.70
CAL.2	5.40	ng/mL	B/T %	1.15
CAL.3	20.00	ng/mL	B/T %	3.77
CAL.4	81.00	ng/mL	B/T %	13.77
CAL.5	210.00	ng/mL	B/T %	30.77

Activité totale / Total activity : 177980 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	4.320 - 6.480	5.41

Date du contrôle
Control Date

11/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

04/07/2022

