

CEA-RIACT

Protocole/Protocol : 1h

Lot N°/Batch # : 479B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-CEA-RIACT	479A	15/02/2021
CT-CEA-RIACT	58A	23/11/2021
DIL-CEA-RIACT	27A	28/10/2021
CAL0-CEA-RIACT	27B	20/03/2021
CAL1-CEA-RIACT	27B	20/03/2021
CAL2-CEA-RIACT	27B	20/03/2021
CAL3-CEA-RIACT	27B	20/03/2021
CAL4-CEA-RIACT	27B	20/03/2021
CAL5-CEA-RIACT	27B	20/03/2021
CONT-CEA-RIACT	27B	20/03/2021
TWEEN-1-3-EL-1FLA	213A	20/03/2023
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 %	0.15
CAL.1	3.10	ng/mL	B/T %	0.81
CAL.2	6.30	ng/mL	B/T %	1.42
CAL.3	20.70	ng/mL	B/T %	3.79
CAL.4	81.00	ng/mL	B/T %	13.79
CAL.5	200.00	ng/mL	B/T %	30.53

Activité totale / Total activity : 188634 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	4.480 - 6.720	5.17

Date du contrôle
Control Date

23/12/2020

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

15/02/2021

