

# CEA-RIACT

Protocole/Protocol : 1h

Lot N°/Batch # : 479A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-CEA-RIACT	479A	15/02/2021
CT-CEA-RIACT	57A	03/08/2021
DIL-CEA-RIACT	27A	28/10/2021
CAL0-CEA-RIACT	27A	20/03/2021
CAL1-CEA-RIACT	27A	20/03/2021
CAL2-CEA-RIACT	27A	20/03/2021
CAL3-CEA-RIACT	27A	20/03/2021
CAL4-CEA-RIACT	27A	20/03/2021
CAL5-CEA-RIACT	27A	20/03/2021
CONT-CEA-RIACT	27A	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 %	0.09
CAL.1	2.90	ng/mL	B/T %	0.79
CAL.2	5.80	ng/mL	B/T %	1.38
CAL.3	19.40	ng/mL	B/T %	4.12
CAL.4	77.00	ng/mL	B/T %	14.28
CAL.5	190.00	ng/mL	B/T %	31.68

Activité totale / Total activity : 184097 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	4.480 - 6.720	4.90

Date du contrôle  
Control Date

22/12/2020

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

15/02/2021

