

# PSA-RIACT

Protocole/Protocol : 2h

Lot N°/Batch # : 458A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-PSA-RIACT	458A	04/07/2022
CT-PSA-RIACT	87A	31/08/2022
CAL0-PSA-RIACT	55A	14/02/2023
CAL1-PSA-RIACT	54A	03/10/2022
CAL2-PSA-RIACT	54A	03/10/2022
CAL3-PSA-RIACT	54A	03/10/2022
CAL4-PSA-RIACT	54A	03/10/2022
CAL5-PSA-RIACT	54A	03/10/2022
CONT-PSA-RIACT	54A	03/10/2022
TWEEN-1-3-EL-1FLA	215A	03/06/2024
PSA-RIA-CT MO	024	

## RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B/T %	0.13
CAL.1	1.10	ng/mL	B/T %	0.78
CAL.2	5.30	ng/mL	B/T %	3.47
CAL.3	20.80	ng/mL	B/T %	13.14
CAL.4	51.00	ng/mL	B/T %	29.28
CAL.5	101.00	ng/mL	B/T %	52.02

Activité totale / Total activity : 196598 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	2.89 - 3.91	3.53

Date du contrôle  
Control Date

13/05/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**04/07/2022**

