

PSA-RIACT

Protocole/Protocol : 2h

Lot N°/Batch # : 451A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-PSA-RIACT	451A	21/12/2021
CT-PSA-RIACT	86A	27/01/2022
CAL0-PSA-RIACT	54A	26/04/2022
CAL1-PSA-RIACT	53B	04/01/2022
CAL2-PSA-RIACT	53B	04/01/2022
CAL3-PSA-RIACT	53B	04/01/2022
CAL4-PSA-RIACT	53A	04/01/2022
CAL5-PSA-RIACT	53A	04/01/2022
CONT-PSA-RIACT	53A	04/01/2022
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 %	0.08
CAL.1	1.00	ng/mL	B/T %	0.76
CAL.2	5.70	ng/mL	B/T %	3.96
CAL.3	21.10	ng/mL	B/T %	14.26
CAL.4	51.00	ng/mL	B/T %	32.33
CAL.5	101.00	ng/mL	B/T %	57.87

Activité totale / Total activity : 186561 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	2.63 - 3.57	3.27

Date du contrôle
Control Date

03/11/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

21/12/2021



DELEUZE Isabelle