

FPSA-RIACT

Protocole/Protocol : 16 - 20h

Lot N°/Batch # : 373A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-FPSA-RIACT	373A	04/07/2022
CT-FPSA-RIACT	69A	14/09/2022
CAL0-FPSA-RIACT	57A	30/12/2022
CAL1-FPSA-RIACT	57B	31/12/2022
CAL2-FPSA-RIACT	57A	30/12/2022
CAL3-FPSA-RIACT	57A	30/12/2022
CAL4-FPSA-RIACT	57A	30/12/2022
CAL5-FPSA-RIACT	57A	30/12/2022
CONT-FPSA-RIACT	57A	30/12/2022
TWEEN-1-3-EL-1FLA	215A	03/06/2024
FPSA-RIA-CT MO	022	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B/T %	0.03
CAL.1	0.48	ng/mL	B/T %	1.55
CAL.2	1.02	ng/mL	B/T %	3.12
CAL.3	2.59	ng/mL	B/T %	7.97
CAL.4	5.20	ng/mL	B/T %	15.29
CAL.5	19.80	ng/mL	B/T %	58.51

Activité totale / Total activity : 174570 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	0.65 - 0.88	0.70

Date du contrôle
Control Date

12/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

04/07/2022

