

IGF1-RIACT

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 233A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-IGF1-RIACT	233A	28/03/2022
CT-IGF1-RIACT	46A	30/11/2022
BUF-IGF1-RIACT	25A	06/04/2022
DIL-IGF1-RIACT	44A	27/06/2022
CAL0-IGF1-RIACT	31A	22/09/2022
CAL1-IGF1-RIACT	31A	22/09/2022
CAL2-IGF1-RIACT	31A	22/09/2022
CAL3-IGF1-RIACT	31A	22/09/2022
CAL4-IGF1-RIACT	31A	22/09/2022
CONT1-IGF1-RIACT	32A	01/11/2022
CONT2-IGF1-RIACT	31B	02/08/2022
TWEEN-1-3-EL-1FLA	215A	03/06/2024
IGF-I-RIACT MO	019	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B/T %	0.26
CAL.1	28.00	ng/mL	B/T %	4.26
CAL.2	141.00	ng/mL	B/T %	17.59
CAL.3	450.00	ng/mL	B/T %	42.13
CAL.4	950.00	ng/mL	B/T %	68.55

Activité totale / Total activity : 189297 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	ng/mL	41.60 - 62.40	47.87
CONTROL 2	ng/mL	280.00 - 420.00	319.14

Date du contrôle
Control Date

05/01/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

28/03/2022

