

# IGF1-RIACT

Protocole/Protocol : 2h + 2h

Lot N°/Batch # : 235A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-IGF1-RIACT	235A	18/07/2022
CT-IGF1-RIACT	46A	30/11/2022
BUF-IGF1-RIACT	26A	02/01/2024
DIL-IGF1-RIACT	45A	23/02/2023
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 %	0.39
CAL.1	28.00	ng/mL	B/T %	4.50
CAL.2	141.00	ng/mL	B/T %	18.24
CAL.3	450.00	ng/mL	B/T %	44.75
CAL.4	950.00	ng/mL	B/T %	69.32

Activité totale / Total activity : 183262 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	ng/mL	41.60 - 62.40	51.48
CONTROL 2	ng/mL	280.00 - 420.00	315.32

Date du contrôle  
Control Date

25/04/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**18/07/2022**

