

# HGH-RIACT

Protocole/Protocol : 2h

Lot N°/Batch # : 271A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-HGH-RIACT	271A	11/03/2022
CT-HGH-RIACT	36A	12/10/2022
CAL0-HGH-RIACT	21A	10/02/2023
CAL1-HGH-RIACT	21A	19/05/2023
CAL2-HGH-RIACT	21A	19/05/2023
CAL3-HGH-RIACT	21A	19/05/2023
CAL4-HGH-RIACT	21A	19/05/2023
CAL5-HGH-RIACT	21A	19/05/2023
CONT-HGH-RIACT	21A	19/05/2023
TWEEN-1-3-EL-1FLA	215A	03/06/2024
HGH-RIA-CT MO	019	

## RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL 0	0.00	ng/mL	B/T %	0.13
CAL 1	0.09	ng/mL	B/T %	0.48
CAL 2	0.93	ng/mL	B/T %	3.54
CAL 3	4.80	ng/mL	B/T %	16.44
CAL 4	12.40	ng/mL	B/T %	36.38
CAL 5	22.50	ng/mL	B/T %	57.09

Activité totale / Total activity : 133445 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	2.720 - 4.080	3.29

Date du contrôle  
Control Date

13/01/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

11/03/2022

