

HGH-RIACT

Protocole/Protocol : 2h

Lot N°/Batch # : 265A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-HGH-RIACT	265A	24/09/2021
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 %	0.07
CAL 1	0.09	ng/mL	B/T %	0.38
CAL 2	0.93	ng/mL	B/T %	3.48
CAL 3	4.80	ng/mL	B/T %	16.20
CAL 4	12.40	ng/mL	B/T %	34.20
CAL 5	22.50	ng/mL	B/T %	58.33

Activité totale / Total activity : 141359 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	2.720 - 3.780	3.27

Date du contrôle
Control Date

26/07/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

24/09/2021

