

OSTEO-US

Protocole/Protocol : 2h

Lot N°/Batch # : 265U

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OSTEO-RIACT	265A	09/12/2021
CT-OSTEO-RIACT	40A	17/01/2022
CAL0-OSTEO-RIACT	20A	31/07/2023
CAL1-OSTEO-RIACT	20A	09/01/2023
CAL2-OSTEO-RIACT	20A	09/01/2023
CAL3-OSTEO-RIACT	20A	09/01/2023
CAL4-OSTEO-RIACT	20A	09/01/2023
CONT-OSTEO-RIACT	20A	09/01/2023
TWEEN-1-3-EL-1FLA	215A	03/06/2024
OSTEO-RIA-CT US MO	015	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B/T %	0.04
CAL.1	4.80	ng/mL	B/T %	0.53
CAL.2	39.00	ng/mL	B/T %	3.79
CAL.3	118.00	ng/mL	B/T %	10.40
CAL.4	280.00	ng/mL	B/T %	25.96

Activité totale / Total activity : 279905 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	14.600 - 22.000	18.63

Date du contrôle
Control Date

13/10/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

09/12/2021

