

# OSTEO-RIACT

Protocole/Protocol : 2h

Lot N°/Batch # : 272A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OSTEO-RIACT	272A	30/06/2022
CT-OSTEO-RIACT	41A	17/10/2022
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 MO	015	

## RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B/T %	0.07
CAL.1	4.80	ng/mL	B/T %	0.54
CAL.2	39.00	ng/mL	B/T %	4.22
CAL.3	118.00	ng/mL	B/T %	12.30
CAL.4	280.00	ng/mL	B/T %	29.03

Activité totale / Total activity : 269907 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	14.600 - 22.000	18.97

Date du contrôle  
Control Date

04/05/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**30/06/2022**

