

# ELSA-PTH

Protocole/Protocol : 16 - 20h

Lot N°/Batch # : 430B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-PTH	430A	02/03/2022
CT-ELSA-PTH	88	29/11/2022
CAL0-ELSA-PTH	34A	07/02/2023
CAL1-ELSA-PTH	33A	14/08/2023
CAL2-ELSA-PTH	33A	14/08/2023
CAL3-ELSA-PTH	33A	14/08/2023
CAL4-ELSA-PTH	33A	14/08/2023
CAL5-ELSA-PTH	33A	14/08/2023
CONT-ELSA-PTH	33A	14/08/2023
TWEEN-1-3-EL-1FLA	215A	03/06/2024
ELSA-PTH MO	025	

## RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	pg/mL	B/T %	0.05
CAL.1	17.00	pg/mL	B/T %	0.60
CAL.2	53.00	pg/mL	B/T %	1.75
CAL.3	160.00	pg/mL	B/T %	5.31
CAL.4	480.00	pg/mL	B/T %	14.80
CAL.5	1 800.00	pg/mL	B/T %	34.07

Activité totale / Total activity : 277742 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	35.200 - 52.800	41.60

Date du contrôle  
Control Date

11/01/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**02/03/2022**

