

ELSA-PTH

Protocole/Protocol : 16 - 20h

Lot N°/Batch # : 430A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-PTH	430A	02/03/2022
CT-ELSA-PTH	87	05/07/2022
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 %	0.06
CAL.1	17.00	pg/mL	B/T %	0.57
CAL.2	53.00	pg/mL	B/T %	1.65
CAL.3	160.00	pg/mL	B/T %	4.95
CAL.4	480.00	pg/mL	B/T %	14.17
CAL.5	1 800.00	pg/mL	B/T %	32.94

Activité totale / Total activity : 273217 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROLA	pg/mL	35.200 - 52.800	43.14

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

