

ELSA-ACTH

Lot N°/Batch # : 433B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-ACTH	433A	09/03/2022
CT-ELSA-ACTH	100	13/12/2022
DIL-ELSA-ACTH	36A	12/10/2022
CAL0-ELSA-ACTH	36A	05/09/2023
CAL1-ELSA-ACTH	35A	24/10/2023
CAL2-ELSA-ACTH	35A	24/10/2023
CAL3-ELSA-ACTH	35A	24/10/2023
CAL4-ELSA-ACTH	35A	24/10/2023
CAL5-ELSA-ACTH	35A	24/10/2023
CONT-ELSA-ACTH	35A	24/10/2023
TWEEN-1-3-EL-1FLA	215A	03/06/2024
ELSA.ACTH 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	19.00	pg/mL	B/T %	0.23
CAL.2	53.00	pg/mL	B/T %	0.62
CAL.3	350.00	pg/mL	B/T %	4.14
CAL.4	880.00	pg/mL	B/T %	9.33
CAL.5	2 000.00	pg/mL	B/T %	22.61

Activité totale / Total activity : 280036 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	35.200 - 52.800	46.89

Date du contrôle
Control Date

10/01/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

09/03/2022

