

ELSA-ACTH

Protocole/Protocol : 18 - 22h

Lot N°/Batch # : 427A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-ACTH	427A	22/09/2021
CT-ELSA-ACTH	99	04/05/2022
DIL-ELSA-ACTH	36A	12/10/2022
CAL0-ELSA-ACTH	35A	23/12/2021
CAL1-ELSA-ACTH	34A	04/02/2022
CAL2-ELSA-ACTH	34A	04/02/2022
CAL3-ELSA-ACTH	34A	04/02/2022
CAL4-ELSA-ACTH	34A	04/02/2022
CAL5-ELSA-ACTH	34A	04/02/2022
CONT-ELSA-ACTH	34A	04/02/2022
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 %	0.03
CAL.1	15.00	pg/mL	B/T %	0.18
CAL.2	49.00	pg/mL	B/T %	0.53
CAL.3	295.00	pg/mL	B/T %	4.04
CAL.4	830.00	pg/mL	B/T %	11.18
CAL.5	2 100.00	pg/mL	B/T %	25.84

Activité totale / Total activity : 293134 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	28.800 - 43.200	39.84

Date du contrôle
Control Date

22/07/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

22/09/2021

