

ELSA-CA125II

Protocole/Protocol : 18 - 22h

Lot N°/Batch # : 436A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-CA125II	436A	24/02/2021
CT-ELSA-CA125II	153	28/04/2021
CAL0-ELSA-CA125II	61A	03/06/2021
CAL1-ELSA-CA125II	67A	02/07/2021
CAL2-ELSA-CA125II	67A	02/07/2021
CAL3-ELSA-CA125II	67A	02/07/2021
CAL4-ELSA-CA125II	67A	02/07/2021
CAL5-ELSA-CA125II	67A	02/07/2021
CONT-ELSA-CA125II	67A	02/07/2021
ELSA-CA-125 2 MO	025	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	U/mL	B/T %	0.06
CAL.1	16.00	U/mL	B/T %	1.02
CAL.2	28.00	U/mL	B/T %	2.03
CAL.3	70.00	U/mL	B/T %	5.35
CAL.4	185.00	U/mL	B/T %	13.49
CAL.5	450.00	U/mL	B/T %	32.32

Activité totale / Total activity : 230743 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	U/mL	88.00 - 132.00	110.73

Date du contrôle
Control Date

28/12/2020

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

24/02/2021

