

ELSA-CA125II

Protocole/Protocol : 18 - 22h

Lot N°/Batch # : 454A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-CA125II	454A	13/07/2022
CT-ELSA-CA125II	156V	13/07/2022
CAL0-ELSA-CA125II	63A	02/12/2022
CAL1-ELSA-CA125II	69A	21/10/2022
CAL2-ELSA-CA125II	69A	21/10/2022
CAL3-ELSA-CA125II	69A	21/10/2022
CAL4-ELSA-CA125II	69A	21/10/2022
CAL5-ELSA-CA125II	69A	21/10/2022
CONT-ELSA-CA125II	69A	21/10/2022
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 %	0.03
CAL.1	15.00	U/mL	B/T %	0.83
CAL.2	28.00	U/mL	B/T %	1.63
CAL.3	74.00	U/mL	B/T %	4.38
CAL.4	185.00	U/mL	B/T %	11.43
CAL.5	500.00	U/mL	B/T %	28.88

Activité totale / Total activity : 234875 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	U/mL	88.00 - 132.00	108.96

Date du contrôle
Control Date

12/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

13/07/2022

