

ELSA-CA19-9

Protocole/Protocol : 3h + 3h

Lot N°/Batch # : 733A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-CA19-9	733A	17/02/2021
CT-ELSA-CA19-9	446	02/06/2021
BUF-ELSA-CA19-9	92A	21/06/2021
DIL-ELSA-CA19-9	218A	10/09/2021
CAL0-ELSA-CA19-9	218A	10/09/2021
CAL1-ELSA-CA19-9	208A	29/09/2021
CAL2-ELSA-CA19-9	208A	29/09/2021
CAL3-ELSA-CA19-9	208A	29/09/2021
CAL4-ELSA-CA19-9	208A	29/09/2021
CAL5-ELSA-CA19-9	208A	29/09/2021
CONT-ELSA-CA19-9	208A	29/09/2021
ELSA CA19-9 MO	039	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	U/mL	B/T %	0.10
CAL.1	13.50	U/mL	B/T %	1.55
CAL.2	26.00	U/mL	B/T %	3.43
CAL.3	54.00	U/mL	B/T %	8.98
CAL.4	125.00	U/mL	B/T %	17.87
CAL.5	270.00	U/mL	B/T %	29.17

Activité totale / Total activity : 184146 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	U/mL	28.00 - 42.00	34.69

Date du contrôle
Control Date

17/12/2020

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

17/02/2021

