

ELSA-CA19-9

Protocole/Protocol : 3h + 16 - 24h

Lot N°/Batch # : 751A ON

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ELSA-CA19-9	751A	06/07/2022
CT-ELSA-CA19-9	460	08/11/2022
BUF-ELSA-CA19-9	94A	26/10/2022
DIL-ELSA-CA19-9	220A	19/01/2023
CAL0-ELSA-CA19-9	220A	19/01/2023
CAL1-ELSA-CA19-9	210B	09/12/2022
CAL2-ELSA-CA19-9	210B	09/12/2022
CAL3-ELSA-CA19-9	210B	09/12/2022
CAL4-ELSA-CA19-9	210B	09/12/2022
CAL5-ELSA-CA19-9	210B	09/12/2022
CONT-ELSA-CA19-9	210B	09/12/2022
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 %	0.23
CAL.1	14.60	U/mL	B/T %	2.35
CAL.2	29.50	U/mL	B/T %	4.84
CAL.3	52.00	U/mL	B/T %	9.11
CAL.4	104.00	U/mL	B/T %	17.40
CAL.5	255.00	U/mL	B/T %	31.71

Activité totale / Total activity : 183356 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	U/mL	30.40 - 45.60	37.70

Date du contrôle
Control Date

04/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

06/07/2022

