

OCFD03-FT4

Protocole/Protocol : 30mn + 1h

Lot N°/Batch # : 376B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCFD03-FT4	376A	27/07/2022
CT-OCFD03-FT4	147B	25/01/2023
BUF-OCFD03-FT4	47A	12/05/2023
CAL0-OCFD03-FT4	18F	03/10/2022
CAL1-OCFD03-FT4	18F	03/10/2022
CAL2-OCFD03-FT4	18F	03/10/2022
CAL3-OCFD03-FT4	18F	03/10/2022
CAL4-OCFD03-FT4	18F	03/10/2022
CAL5-OCFD03-FT4	18F	03/10/2022
CAL6-OCFD03-FT4	18F	03/10/2022
CONT-OCFD03-FT4	18F	03/10/2022
RG-FT4 MO	024	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	pg/mL	B0/T %	56.26
CAL.1	2.40	pg/mL	B/B0	81.89
CAL.2	4.80	pg/mL	B/B0	67.80
CAL.3	12.00	pg/mL	B/B0	57.56
CAL.4	23.70	pg/mL	B/B0	42.70
CAL.5	44.00	pg/mL	B/B0	31.40
CAL.6	66.00	pg/mL	B/B0	25.44

Activité totale / Total activity : 74861 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	10.40 - 15.60	13.59

Date du contrôle
Control Date

17/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

27/07/2022

