

OCFD03-FT4-K

Protocole/Protocol : 45mn

Lot N°/Batch # : 368A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCFD03-FT4	368A	15/12/2021
CT-OCFD03-FT4	144A	03/08/2022
CAL0 OCFD03-FT4-K	20E	24/01/2022
CAL1 OCFD03-FT4-K	20E	24/01/2022
CAL2 OCFD03-FT4-K	20E	24/01/2022
CAL3 OCFD03-FT4-K	20E	24/01/2022
CAL4 OCFD03-FT4-K	20E	24/01/2022
CAL5 OCFD03-FT4-K	20E	24/01/2022
CAL6 OCFD03-FT4-K	20E	24/01/2022
CONT OCFD03-FT4-K	20E	24/01/2022
OCFD03-FT4-K MO	002	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	pg/mL	B0/T %	52.74
CAL.1	3.00	pg/mL	B/B0	73.06
CAL.2	6.00	pg/mL	B/B0	60.71
CAL.3	10.50	pg/mL	B/B0	48.35
CAL.4	20.00	pg/mL	B/B0	32.80
CAL.5	42.00	pg/mL	B/B0	17.43
CAL.6	68.00	pg/mL	B/B0	9.06

Activité totale / Total activity : 75452 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	7.36 - 11.04	9.12

Date du contrôle
Control Date

06/10/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

15/12/2021

