

# OCFD03-FT4

Protocole/Protocol : 30mn + 1h

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
BUF-OCFD03-FT4	47A	12/05/2023
CAL0-OCFD03-FT4	18E	24/01/2022
CAL1-OCFD03-FT4	18E	24/01/2022
CAL2-OCFD03-FT4	18E	24/01/2022
CAL3-OCFD03-FT4	18E	24/01/2022
CAL4-OCFD03-FT4	18E	24/01/2022
CAL5-OCFD03-FT4	18E	24/01/2022
CAL6-OCFD03-FT4	18E	24/01/2022
CONT-OCFD03-FT4	18E	24/01/2022
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 %	55.42
CAL.1	2.40	pg/mL	B/B0	84.67
CAL.2	4.80	pg/mL	B/B0	71.83
CAL.3	12.00	pg/mL	B/B0	55.49
CAL.4	23.70	pg/mL	B/B0	42.10
CAL.5	44.00	pg/mL	B/B0	30.46
CAL.6	66.00	pg/mL	B/B0	22.38

Activité totale / Total activity : 75331 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	10.40 - 15.60	14.91

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

