

OCFD03-FT4

Protocole/Protocol : 30mn + 1h

Lot N°/Batch # : 369A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCFD03-FT4	369A	12/01/2022
CT-OCFD03-FT4	144B	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 %	55.45
CAL.1	2.40	pg/mL	B/B0	80.90
CAL.2	4.80	pg/mL	B/B0	69.85
CAL.3	12.00	pg/mL	B/B0	55.27
CAL.4	23.70	pg/mL	B/B0	39.58
CAL.5	44.00	pg/mL	B/B0	29.93
CAL.6	66.00	pg/mL	B/B0	23.40

Activité totale / Total activity : 74464 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	10.400 - 15.600	14.40

Date du contrôle
Control Date

03/11/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

12/01/2022



DELEUZE Isabelle