

OCFH07-FT3

Protocole/Protocol : 2h

Lot N°/Batch # : 368A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCFH07-FT3	368A	15/12/2021
CT-OCFH07-FT3	109A	02/03/2022
CAL0-OCFH07-FT3	22C	03/08/2022
CAL1-OCFH07-FT3	22C	03/08/2022
CAL2-OCFH07-FT3	22C	03/08/2022
CAL3-OCFH07-FT3	22C	03/08/2022
CAL4-OCFH07-FT3	22C	03/08/2022
CAL5-OCFH07-FT3	22C	03/08/2022
CAL6-OCFH07-FT3	22C	03/08/2022
CONT-OCFH07-FT3	22C	03/08/2022
RG-FT3 MO	024	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	pg/mL	B0/T %	46.64
CAL.1	0.86	pg/mL	B/B0	85.83
CAL.2	2.03	pg/mL	B/B0	74.40
CAL.3	4.10	pg/mL	B/B0	58.04
CAL.4	9.50	pg/mL	B/B0	39.31
CAL.5	25.90	pg/mL	B/B0	18.58
CAL.6	52.00	pg/mL	B/B0	7.58

Activité totale / Total activity : 68653 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/ml	3.120 - 4.680	3.64

Date du contrôle
Control Date

07/10/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

15/12/2021

