

OCFH07-FT3

Protocole/Protocol : 2h

Lot N°/Batch # : 376A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCFH07-FT3	376A	27/07/2022
CT-OCFH07-FT3	111A	03/08/2022
CAL0-OCFH07-FT3	22E	19/04/2023
CAL1-OCFH07-FT3	22E	19/04/2023
CAL2-OCFH07-FT3	22E	19/04/2023
CAL3-OCFH07-FT3	22E	19/04/2023
CAL4-OCFH07-FT3	22E	19/04/2023
CAL5-OCFH07-FT3	22E	19/04/2023
CAL6-OCFH07-FT3	22E	19/04/2023
CONT-OCFH07-FT3	22E	19/04/2023
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 %	43.52
CAL.1	0.86	pg/mL	B/B0	85.21
CAL.2	2.03	pg/mL	B/B0	73.72
CAL.3	4.10	pg/mL	B/B0	56.78
CAL.4	9.50	pg/mL	B/B0	37.93
CAL.5	25.90	pg/mL	B/B0	18.10
CAL.6	52.00	pg/mL	B/B0	7.42

Activité totale / Total activity : 66490 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/ml	3.12 - 4.68	3.73

Date du contrôle
Control Date

19/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

27/07/2022

