

# OCFD03-FT4

Protocole/Protocol : 30mn + 1h

Lot N°/Batch # : 376A

| Réactifs<br>Reagents | Lot N°<br>Batch # | Péremption<br>Expiry Date |
|----------------------|-------------------|---------------------------|
| TRA-OCFD03-FT4       | 376A              | 27/07/2022                |
| CT-OCFD03-FT4        | 147A              | 25/01/2023                |
| BUF-OCFD03-FT4       | 47A               | 12/05/2023                |
| CAL0-OCFD03-FT4      | 18F               | 03/10/2022                |
| CAL1-OCFD03-FT4      | 18F               | 03/10/2022                |
| CAL2-OCFD03-FT4      | 18F               | 03/10/2022                |
| CAL3-OCFD03-FT4      | 18F               | 03/10/2022                |
| CAL4-OCFD03-FT4      | 18F               | 03/10/2022                |
| CAL5-OCFD03-FT4      | 18F               | 03/10/2022                |
| CAL6-OCFD03-FT4      | 18F               | 03/10/2022                |
| CONT-OCFD03-FT4      | 18F               | 03/10/2022                |
| RG-FT4 MO            | 024               |                           |

## RESULTATS / RESULTS

| Paramètre<br>Parameter | Conc.<br>Conc. | Unité<br>Unit | Réponse<br>Response | Valeur mesurée<br>Measured value |
|------------------------|----------------|---------------|---------------------|----------------------------------|
| CAL.0                  | 0.00           | pg/mL         | B0/T %              | 53.90                            |
| CAL.1                  | 2.40           | pg/mL         | B/B0                | 86.20                            |
| CAL.2                  | 4.80           | pg/mL         | B/B0                | 73.72                            |
| CAL.3                  | 12.00          | pg/mL         | B/B0                | 61.53                            |
| CAL.4                  | 23.70          | pg/mL         | B/B0                | 45.36                            |
| CAL.5                  | 44.00          | pg/mL         | B/B0                | 33.90                            |
| CAL.6                  | 66.00          | pg/mL         | B/B0                | 25.99                            |

Activité totale / Total activity : 75746 cpm

| Contrôles<br>Controls | Unité<br>Unit | Valeurs attendues<br>Expected values | Valeur mesurée<br>Measured value |
|-----------------------|---------------|--------------------------------------|----------------------------------|
| CONTROL               | pg/mL         | 10.40 - 15.60                        | 14.77                            |

Date du contrôle  
Control Date

18/05/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**27/07/2022**

