

# OCPG07-T4

Protocole/Protocol : 45mn

Lot N°/Batch # : 283A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCPG07-T4	283A	02/11/2021
CT-OCPG07-T4	76A	30/03/2022
CAL0-OCPG07-T4	18D	22/04/2022
CAL1-OCPG07-T4	18D	22/04/2022
CAL2-OCPG07-T4	18D	22/04/2022
CAL3-OCPG07-T4	18D	22/04/2022
CAL4-OCPG07-T4	18D	22/04/2022
CAL5-OCPG07-T4	18D	22/04/2022
CONT-OCPG07-T4	18D	22/04/2022
RG-T4 MO	021	

## RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B0/T %	65.83
CAL.1	22.60	ng/mL	B/B0	49.47
CAL.2	49.00	ng/mL	B/B0	35.15
CAL.3	100.00	ng/mL	B/B0	20.25
CAL.4	196.00	ng/mL	B/B0	12.75
CAL.5	340.00	ng/mL	B/B0	8.50

Activité totale / Total activity : 72220 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	64.000 - 96.000	76.23

Date du contrôle  
Control Date

19/08/2021

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**02/11/2021**

