

# T4-US

Protocole/Protocol : 45mn

Lot N°/Batch # : 292U

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-T4-US	292U	
CT-T4-US	77U	
CAL0-T4-US	18X	
CAL1-T4-US	18X	
CAL2-T4-US	18X	
CAL3-T4-US	18X	
CAL4-T4-US	18X	
CAL5-T4-US	18X	
CT-T4-US	18X	
RG-T4-US MO	004	

## RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B0/T %	77.38
CAL.1	22.60	ng/mL	B/B0	53.67
CAL.2	49.00	ng/mL	B/B0	34.63
CAL.3	100.00	ng/mL	B/B0	20.80
CAL.4	196.00	ng/mL	B/B0	13.15
CAL.5	340.00	ng/mL	B/B0	8.65

Activité totale / Total activity : 67810 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	ng/mL	64.00 - 96.00	84.50

Date du contrôle  
Control Date

27/04/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

12/07/2022

