

OCPM03-HCG

Protocole/Protocol : 30mn + 1h

Lot N°/Batch # : 237A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-OCPM03-HCG	237A	12/11/2021
CT-OCPM03-HCG	56A	17/02/2022
BUF-OCPM03-HCG	17A	16/08/2023
CAL0-OCPM03-HCG	23B	16/06/2022
CAL1-OCPM03-HCG	23B	16/06/2022
CAL2-OCPM03-HCG	23B	16/06/2022
CAL3-OCPM03-HCG	23B	16/06/2022
CAL4-OCPM03-HCG	23B	16/06/2022
CAL5-OCPM03-HCG	23B	16/06/2022
CAL6-OCPM03-HCG	23B	16/06/2022
CAL7-OCPM03-HCG	23B	16/06/2022
CONT-OCPM03-HCG	23B	16/06/2022
BLISTER-5-WASH	25B	28/02/2022
RG-HCG MO	017	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	mUI/mL	B/T %	0.08
CAL.1	4.70	mUI/mL	B/T %	0.49
CAL.2	10.80	mUI/mL	B/T %	1.08
CAL.3	31.00	mUI/mL	B/T %	2.82
CAL.4	114.00	mUI/mL	B/T %	10.17
CAL.5	300.00	mUI/mL	B/T %	25.24
CAL.6	640.00	mUI/mL	B/T %	43.33
CAL.7	1 080.00	mUI/mL	B/T %	52.21

Activité totale / Total activity : 151106 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	mUI/mL	96.00 - 144.00	122.45

Date du contrôle
Control Date

20/08/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

12/11/2021

