

R-GM-100

Protocole/Protocol : 1h

Lot N°/Batch # : 59A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-ANDROSTENEDIONE	59A	30/09/2021
CT-ANDROSTENEDIONE	10A	24/02/2022
CAL0-ANDROSTENEDIONE	17A	26/10/2022
CAL1-ANDROSTENEDIONE	17A	26/10/2022
CAL2-ANDROSTENEDIONE	17A	26/10/2022
CAL3-ANDROSTENEDIONE	17A	26/10/2022
CAL4-ANDROSTENEDIONE	17A	26/10/2022
CAL5-ANDROSTENEDIONE	18A	26/10/2022
CONT1-ANDROSTENEDION	17A	26/10/2022
CONT2-ANDROSTENEDION	17A	26/10/2022
TWEEN-1-3-EL-1FLA	214A	18/03/2024
ANDROSTENEDIONE MO	003	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	ng/mL	B0/T %	27.73
CAL.1	0.10	ng/mL	B/B0	79.61
CAL.2	0.32	ng/mL	B/B0	63.88
CAL.3	0.92	ng/mL	B/B0	43.61
CAL.4	2.70	ng/mL	B/B0	25.17
CAL.5	11.50	ng/mL	B/B0	9.72

Activité totale / Total activity : 52424 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL 1	ng/mL	0.610 - 0.910	0.71
CONTROL 2	ng/mL	3.140 - 4.700	4.06

Date du contrôle
Control Date

19/07/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

30/09/2021

