

IRMA-HCT

Protocole/Protocol : 20 - 24h

Lot N°/Batch # : 499A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-IRMA-HCT	499A	06/09/2021
CT-IRMA-HCT	40A	14/09/2021
CAL0-IRMA-HCT	20A	04/04/2022
CAL1-IRMA-HCT	27A	05/10/2021
CAL2-IRMA-HCT	27A	05/10/2021
CAL3-IRMA-HCT	27A	05/10/2021
CAL4-IRMA-HCT	27A	05/10/2021
CAL5-IRMA-HCT	27A	05/10/2021
CONT-IRMA-HCT	27A	05/10/2021
BLISTER-5-WASH	25B	28/02/2022
IRMA-HCT MO	020	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	pg/mL	B/T %	0.01
CAL.1	7.90	pg/mL	B/T %	0.16
CAL.2	31.00	pg/mL	B/T %	0.56
CAL.3	110.00	pg/mL	B/T %	2.79
CAL.4	710.00	pg/mL	B/T %	21.88
CAL.5	1 400.00	pg/mL	B/T %	42.60

Activité totale / Total activity : 272841 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	9.280 - 13.920	11.66
AS014	pg/mL	4.80 - 9.00	7.07

Date du contrôle
Control Date

27/07/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

06/09/2021

