

# IRMA-HCT

Protocole/Protocol : 20 - 24h

Lot N°/Batch # : 511A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-IRMA-HCT	511A	21/02/2022
CT-IRMA-HCT	41A	04/07/2022
CAL0-IRMA-HCT	20A	04/04/2022
CAL1-IRMA-HCT	28A	13/07/2023
CAL2-IRMA-HCT	28A	13/07/2023
CAL3-IRMA-HCT	28A	13/07/2023
CAL4-IRMA-HCT	28A	13/07/2023
CAL5-IRMA-HCT	28A	13/07/2023
CONT-IRMA-HCT	28A	13/07/2023
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 %	0.02
CAL.1	6.20	pg/mL	B/T %	0.17
CAL.2	25.00	pg/mL	B/T %	0.60
CAL.3	105.00	pg/mL	B/T %	3.09
CAL.4	650.00	pg/mL	B/T %	22.42
CAL.5	1 200.00	pg/mL	B/T %	42.02

Activité totale / Total activity : 273458 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	7.04 - 10.56	9.53
AS014	pg/mL	4.80 - 9.00	7.04

Date du contrôle  
Control Date

10/01/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

21/02/2022

