

# IRMA-HCT

Protocole/Protocol : 20 - 24h

Lot N°/Batch # : 520A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-IRMA-HCT	520A	30/06/2022
CT-IRMA-HCT	41A	04/07/2022
CAL0-IRMA-HCT	21V	30/06/2022
CAL1-IRMA-HCT	29A	13/07/2023
CAL2-IRMA-HCT	29A	13/07/2023
CAL3-IRMA-HCT	29A	13/07/2023
CAL4-IRMA-HCT	29A	13/07/2023
CAL5-IRMA-HCT	29A	13/07/2023
CONT-IRMA-HCT	29A	13/07/2023
BLISTER-3-WASH	26A	31/03/2025
IRMA-HCT MO	021	

## 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	7.40	pg/mL	B/T %	0.19
CAL.2	27.50	pg/mL	B/T %	0.69
CAL.3	113.00	pg/mL	B/T %	3.45
CAL.4	680.00	pg/mL	B/T %	23.49
CAL.5	1 230.00	pg/mL	B/T %	44.97

Activité totale / Total activity : 280315 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	pg/mL	8.80 - 13.20	10.86
AS014	pg/mL	4.80 - 9.00	6.98

Date du contrôle  
Control Date

17/05/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

**30/06/2022**

