

BI-INS-IRMA

Protocole/Protocol : 2h

Lot N°/Batch # : 266A

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
TRA-BI-INS-IRMA	266A	04/07/2022
CT-BI-INS-IRMA	47A	10/10/2023
RIP-BI-INS-IRMA	16A	04/03/2024
CAL0-BI-INS-IRMA	23A	16/02/2023
CAL1-BI-INS-IRMA	24A	01/06/2023
CAL2-BI-INS-IRMA	24A	01/06/2023
CAL3-BI-INS-IRMA	24A	01/06/2023
CAL4-BI-INS-IRMA	24A	01/06/2023
CAL5-BI-INS-IRMA	24A	01/06/2023
CAL6-BI-INS-IRMA	24A	01/06/2023
CONT-BI-INS-IRMA	16A	02/05/2023
SOL-WASH-R3	16A	23/09/2022
BI-INSULIN IRMA MO	019	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL.0	0.00	µUI/mL	cpm	326.24
CAL.1	1.00	µUI/mL	B-B0/T	0.11
CAL.2	5.00	µUI/mL	B-B0/T	0.87
CAL.3	19.00	µUI/mL	B-B0/T	4.08
CAL.4	100.00	µUI/mL	B-B0/T	18.95
CAL.5	250.00	µUI/mL	B-B0/T	41.46
CAL.6	450.00	µUI/mL	B-B0/T	63.65

Activité totale / Total activity : 245091 cpm

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CONTROL	µUI/mL	17.00 - 25.60	21.17

Date du contrôle
Control Date

11/05/2022

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

04/07/2022

