

P3NP-ELISA

Protocole/Protocol : 3h

Lot N°/Batch # : 13B

Réactifs Reagents	Lot N° Batch #	Péremption Expiry Date
CONJ N Terminal Pro	13A	27/04/2023
MICROPLATE N Termina	12A	06/02/2023
CAL0-DIL N Terminal	10A	07/02/2023
CAL1 N Terminal Pro	11A	28/03/2023
CAL2 N Terminal Pro	11A	28/03/2023
CAL3 N Terminal Pro	11A	28/03/2023
CAL4 N Terminal Pro	11A	28/03/2023
CAL5 N Terminal Pro	11A	28/03/2023
CONT1 N Terminal Pro	11A	28/03/2023
CONT2 N Terminal Pro	11A	28/03/2023
BLISTER-3-WASH	26A	31/03/2025
TWEEN-1-3-EL-1FLA	215A	03/06/2024
SUBS-TMB	55A	08/07/2023
STOP-SOL	29A	01/11/2022
N Terminal Procollag	007	

RESULTATS / RESULTS

Paramètre Parameter	Conc. Conc.	Unité Unit	Réponse Response	Valeur mesurée Measured value
CAL 0	0.00	µg/L	DO	0.06
CAL 1	3.00	µg/L	DO	0.40
CAL 2	6.30	µg/L	DO	0.72
CAL 3	12.00	µg/L	DO	1.24
CAL 4	24.00	µg/L	DO	1.91
CAL 5	35.00	µg/L	DO	2.30

Contrôles Controls	Unité Unit	Valeurs attendues Expected values	Valeur mesurée Measured value
CRT1	µg/L	6.0 - 9.0	7.55
CRT2	µg/L	15.9 - 23.9	20.34

Date du contrôle
Control Date

11/05/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

01/11/2022

