

P3NP-EL-US

Protocole/Protocol : 3h

Lot N°/Batch # : 14V

| Réactifs Reagents | Lot N° Batch # |
|----------------------|-------------------|
| CONJ N Terminal Proc | 14V |
| MICROPLATE N Termina | 13V |
| CAL0-DIL N Terminal | 11V |
| CAL1 N Terminal Proc | 12V |
| CAL2 N Terminal Proc | 12V |
| CAL3 N Terminal Proc | 12V |
| CAL4 N Terminal Proc | 12V |
| CAL5 N Terminal Proc | 12V |
| CONT1 N Terminal Pro | 12V |
| CONT2 N Terminal Pro | 12V |
| TWEEN-20-US | 213V |
| BLISTER-3-WASH | 26A |
| SUBS-TMB-US | 55V |
| STOP-SOL-US | 30V |
| N Terminal Procollag | 009 |

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 | 2.97 | µg/L | DO | 0.46 |
| CAL 2 | 6.20 | µg/L | DO | 0.88 |
| CAL 3 | 12.50 | µg/L | DO | 1.52 |
| CAL 4 | 24.50 | µg/L | DO | 2.24 |
| CAL 5 | 33.00 | µg/L | DO | 2.59 |

| Contrôles Controls | Unité Unit | Valeurs attendues Expected values | Valeur mesurée Measured value |
|-----------------------|---------------|--------------------------------------|----------------------------------|
| CRT1 | µg/L | 5.92 - 8.88 | 8.16 |
| CRT2 | µg/L | 16.000 - 24.000 | 20.98 |

Date du contrôle
Control Date

06/09/2021

Visa du responsable du contrôle qualité
Quality control manager visa

Date de péremption
Expiry date

20/03/2023

