

# P3NP-EL-US

Protocole/Protocol : 3h

Lot N°/Batch # : 15V

| Réactifs<br>Reagents | Lot N°<br>Batch # |
|----------------------|-------------------|
| CONJ-P3NP-EL-US      | 15V               |
| MICROPLATE-P3NP-EL-U | 15V               |
| DIL CAL0-P3NP-EL-US  | 15V               |
| CAL1-P3NP-EL-US      | 15V               |
| CAL2-P3NP-EL-US      | 15V               |
| CAL3-P3NP-EL-US      | 15V               |
| CAL4-P3NP-EL-US      | 15V               |
| CAL5-P3NP-EL-US      | 15V               |
| CONT1-P3NP-EL-US     | 15V               |
| CONT2-P3NP-EL-US     | 15V               |
| TWEEN 20             | 215VP             |
| BLISTER-3-WASH       | 26A               |
| SUBS TMB             | 57VP              |
| STOP SOLN            | 29VP              |
| P3NP US MO           | 010               |

## RESULTATS / RESULTS

| Paramètre<br>Parameter | Conc.<br>Conc. | Unité<br>Unit | Réponse<br>Response | Valeur mesurée<br>Measured value |
|------------------------|----------------|---------------|---------------------|----------------------------------|
| CAL 0                  | 0.00           | µg/L          | DO                  | 0.06                             |
| CAL 1                  | 2.95           | µg/L          | DO                  | 0.45                             |
| CAL 2                  | 6.10           | µg/L          | DO                  | 0.82                             |
| CAL 3                  | 12.10          | µg/L          | DO                  | 1.41                             |
| CAL 4                  | 22.60          | µg/L          | DO                  | 2.07                             |
| CAL 5                  | 32.00          | µg/L          | DO                  | 2.53                             |

| Contrôles<br>Controls | Unité<br>Unit | Valeurs attendues<br>Expected values | Valeur mesurée<br>Measured value |
|-----------------------|---------------|--------------------------------------|----------------------------------|
| CRT1                  | µg/L          | 5.84 - 8.76                          | 7.63                             |
| CRT2                  | µg/L          | 15.4 - 23.2                          | 19.40                            |

Date du contrôle  
Control Date

03/08/2022

Visa du responsable du contrôle qualité  
Quality control manager visa

Date de péremption  
Expiry date

30/04/2024

