

|                 |             |
|-----------------|-------------|
| Product Number: | 101477      |
| Lot Number:     | 504658      |
| Expiration:     | 24-Mar-2026 |
| Storage:        | 2-8°C       |

| Component                     | Part Number | Lot Number | Control Lot Number |
|-------------------------------|-------------|------------|--------------------|
| Sample Diluent A              | 101472      | 501527     | 501527A            |
| Sample Diluent B              | 101473      | 426818     | N/A                |
| Sample Diluent C              | 101474      | 511201     | N/A                |
| Sample Diluent D              | 101475      | 510607     | N/A                |
| Sample Diluent E <sup>1</sup> | 101476      | 516704     | N/A                |

Laboratory Analysis

| Characteristic | Result | pH Target   |
|----------------|--------|-------------|
| pH Measurement | Pass   | 7.35 - 7.45 |

| Characteristic | Result | pH Target                |
|----------------|--------|--------------------------|
| pH Measurement | Pass   | 8.95 - 9.05 <sup>1</sup> |

<sup>1</sup> Sample Diluent E (101476) pH target differs from the all other components  
Note: Result corresponds to each component