

|                 |             |
|-----------------|-------------|
| Product Number: | 101477      |
| Lot Number:     | 504727      |
| Expiration:     | 15-Jul-2026 |
| Storage:        | 2-8°C       |

| Component                     | Part Number | Lot Number | Control Lot Number |
|-------------------------------|-------------|------------|--------------------|
| Sample Diluent A              | 101472      | 501527     | 501527A            |
| Sample Diluent B              | 101473      | 521616     | N/A                |
| Sample Diluent C              | 101474      | 521712     | N/A                |
| Sample Diluent D              | 101475      | 521907     | N/A                |
| Sample Diluent E <sup>1</sup> | 101476      | 525108     | 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

Review/Approval

Dhwani Patel

QA Specialist II, QA

 18 Sep 2025

Name

Title

Signature/Date