

|                 |             |
|-----------------|-------------|
| Product Number: | 101477      |
| Lot Number:     | 504033      |
| Expiration:     | 23-Mar-2024 |
| Storage:        | 2-8°C       |

| Component                     | Part Number | Lot Number |
|-------------------------------|-------------|------------|
| Sample Diluent A              | 101472      | 329408     |
| Sample Diluent B              | 101473      | 308223     |
| Sample Diluent C              | 101474      | 328202     |
| Sample Diluent D              | 101475      | 326107     |
| Sample Diluent E <sup>1</sup> | 101476      | 328612     |

**Laboratory Analysis**

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

*Sonia Khillar*

*Senior QC Analyst*

*Sonia Khillar* 31 OCT 2023

Name

Title

Signature/Date