

|                 |            |
|-----------------|------------|
| Product Number: | 101353     |
| Lot Number:     | 300909     |
| Expiration:     | 9-Jan-2024 |
| Storage:        | 2 - 8°C    |

## Laboratory Analysis

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

Review/Approval

|               |            |                        |
|---------------|------------|------------------------|
| Sonia Khillan | Ac Manager | Sonia/20/1/17 JAN 2023 |
| Name          | Title      | Signature/Date         |