

Product Number:	101476
Lot Number:	201811
Expiration:	18-Jan-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95-9.05

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 28 Mar 2027
Signature/Date