

Product Number:	101476
Lot Number:	305404
Expiration:	23-Feb-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95 - 9.05

Review/Approval

Sonia Khillsn
Name

QC Manager
Title

Sonia Khillsn / 09MAR2023
Signature/Date