

Product Number:	101954
Lot Number:	306808
Expiration:	9-Mar-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 21 MAR 2023
Signature/Date