

Product Number:	101476
Lot Number:	322906
Expiration:	17-Aug-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.96	8.95 - 9.05	PASS

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 01SEP2023
Signature/Date