

Product Number:	103360
Lot Number:	327911
Expiration:	6-Oct-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.98	7.95 - 8.05	PASS

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 11 OCT 2023
Signature/Date