

Product Number:	101956
Lot Number:	327509
Expiration:	2-Oct-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.37	7.35 - 7.45	PASS

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 11 OCT 2023

Signature/Date