

Product Number:	101956
Lot Number:	310016
Expiration:	10-Apr-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 11MAY2023
Signature/Date