

Product Number:	101356
Lot Number:	234208
Expiration:	8-Dec-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan QC Manager Sonia Khillan / 15DEC2022

Name

Title

Signature/Date