

Product Number:	101357
Lot Number:	233907
Expiration:	5-Dec-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 12/16/2022

Signature/Date