

Product Number:	100457
Lot Number:	227911
Expiration:	6-Oct-2023
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 11/3/2022
Signature/Date