

Product Number:	100457
Lot Number:	320106
Expiration:	20-Jul-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 01/AUG/2023
Signature/Date