

Product Number:	101359
Lot Number:	309413
Expiration:	4-Apr-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

<u>Sonia Khillan</u>	<u>QC Manager</u>	<u>Sonia Khillan 10MAY2023</u>
Name	Title	Signature/Date