

Product Number:	100458
Lot Number:	306809
Expiration:	9-Mar-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

<i>Sonia Khillan</i>	<i>QC Manager</i>	<i>Sonia Khillan</i> 21MAR2023
Name	Title	Signature/Date