

Product Number:	100453
Lot Number:	217505
Expiration:	24-Jun-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval

*Sonia Khilkin*

Name

QC

Title

*Sonia Kella / 09 AUG 2022*

Signature/Date