

Product Number:	100453
Lot Number:	306010
Expiration:	1-Mar-2024
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan 07 APR 2023*  
Signature/Date