

Product Number:	100453
Lot Number:	134612
Expiration:	12-Dec-2022
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 20 Apr 2022  
Signature/Date