

Product Number:	100457
Lot Number:	210316
Expiration:	13-Apr-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 20Apr2022  
Signature/Date