

Product Number:	101362
Lot Number:	208307
Expiration:	24-Mar-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 28 Mar 2022  
Signature/Date