

Product Number:	100458
Lot Number:	210310
Expiration:	13-Apr-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 20 Apr 2022  
Signature/Date