

Product Number:	101355
Lot Number:	206115
Expiration:	2-Mar-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Richa Shrestha

QC Supervisor

Name

Title

 29Apr2022  
Signature/Date