

Product Number:	101356
Lot Number:	206118
Expiration:	2-Mar-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 08 Mar 2022

Name

Title

Signature/Date