

Product Number:	100376
Lot Number:	208905
Expiration:	30-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 01Apr 2022

Name

Title

Signature/Date