

Product Number:	103361
Lot Number:	210908
Expiration:	19-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 20Apr 2022
Signature/Date