

Product Number:	103358
Lot Number:	032603
Expiration:	21-Nov-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.9-8.1

Review/Approval

Shivani Goel
Name

Manager, QC
Title

Shivani Goel 01 Dec 2020
Signature/Date