

Product Number:	101472
Lot Number:	024418
Expiration:	31-Aug-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 01 Dec 2020
Signature/Date