

Product Number:	101476
Lot Number:	029511
Expiration:	21-Oct-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 22NOV
Signature/Date 2020