

Product Number:	102913
Lot Number:	124610
Expiration:	3-Sep-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.90-8.10

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 15 Sep 2021
Signature/Date