

Product Number:	103359
Lot Number:	126610
Expiration:	23-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 27sep2021
Signature/Date