

Product Number:	101631
Lot Number:	126908
Expiration:	26-Sep-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.30-7.50

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel (20 Oct 2021)

Signature/Date