

Product Number:	101473
Lot Number:	115227
Expiration:	1-Jun-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 04 Aug 2024

Signature/Date