

Product Number:	101474
Lot Number:	133904
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Gael

Name

QC

Title

Shivani Gael 14 Dec 2021

Signature/Date