

Product Number:	102259
Lot Number:	129305
Expiration:	20-Oct-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 27Oct2021

Signature/Date