

Product Number:	101472
Lot Number:	115921
Expiration:	8-Jun-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shwani Goel

Name

QC Supervisor

Title

Shwani Goel 15 Jun 2022

Signature/Date