

Product Number:	101359
Lot Number:	120409
Expiration:	23-Jul-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 03Aug
Signature/Date 2021