

Product Number:	101359
Lot Number:	134925
Expiration:	15-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivam Goel
Name

QC

Title

Shivam Goel 21Jan2022
Signature/Date