

Product Number:	101357
Lot Number:	134611
Expiration:	12-Dec-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 08 Mar 2022

Name

Title

Signature/Date