

Product Number:	101954
Lot Number:	035808
Expiration:	23-Dec-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Manager, QC

Shivani Goel 15 Jan 2021

Name

Title

Signature/Date