

Product Number:	101954
Lot Number:	122709
Expiration:	15-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

QC

Shivani Goel 21 Sep 2021

Name

Title

Signature/Date