

Product Number:	101362
Lot Number:	031129
Expiration:	6-Nov-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Manager, QC

Shivani Goel 23 Nov 2020

Name

Title

Signature/Date