

Product Number:	101352
Lot Number:	031130
Expiration:	6-Nov-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval Shivani Goel QC Manager Shivani Goel 24 Mar 2021
Name Title Signature/Date