

Product Number:	101362
Lot Number:	105402
Expiration:	23-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval Shivani Goel QC Manager Shivani Goel 04 Mar 2021
Name Title Signature/Date