

Product Number:	101357
Lot Number:	100710
Expiration:	7-Jan-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval Shivani Goel QC Manager Shivani Goel 30 Mar 2021
Name Title Signature/Date