

Product Number:	100453
Lot Number:	120810
Expiration:	26-Jul-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval Shivani Goel QC Shivani Goel 10 Aug 2021
Name Title Signature/Date