

Product Number:	101956
Lot Number:	121717
Expiration:	5-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 26 Aug 2021
Signature/Date