

Product Number:	101476
Lot Number:	019026
Expiration:	8-Jul-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Manager, QC

Name

Title

Shivani Goel 20 Oct 2020

Signature/Date