

Product Number:	102002
Lot Number:	019801
Expiration:	16-Jul-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

Manager, QC
Title

Shivani Goel 05Feb2021
Signature/Date