

Product Number:	101353
Lot Number:	119417
Expiration:	13-Jul-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

Quality Control
Title

Shivani Goel 04 Aug 2021
Signature/Date