

Product Number:	101353
Lot Number:	111312
Expiration:	23-Apr-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

Quality Control
Title

Shivani Goel 03 May 2021
Signature/Date