

Product Number:	101353
Lot Number:	117117
Expiration:	20-Jul-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

Quality Control
Title

Shivani Goel 08 Jul 2022
Signature/Date