

Product Number:	101359
Lot Number:	106318
Expiration:	4-Mar-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

Quality Control
Title

Shivani Goel 24 Mar 2021
Signature/Date