

Product Number:	101359
Lot Number:	117115
Expiration:	20-Jun-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 25-Jun-2021
Signature/Date