

Product Number:	103361
Lot Number:	106704
Expiration:	8-Mar-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH	Pass

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 24 Mar 2021
Signature/Date