

Product Number:	102914
Lot Number:	127502
Expiration:	2-Oct-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH	Pass

Review/Approval

Shivani Goel

Name

QC

Shivani Goel 06Oct2021

Title

Signature/Date