

Product Number:	103646
Lot Number:	124611
Expiration:	3-Sep-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 07 Sep 2021

Signature/Date