

Product Number:	103516
Lot Number:	113505
Expiration:	15-May-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

QC

Shivani Goel 01 Dec 2021

Name

Title

Signature/Date