

Product Number:	101362
Lot Number:	128806
Expiration:	15-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 26 Oct 2021

Signature/Date