

Product Number:	101678
Lot Number:	128802
Expiration:	15-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 19 Oct 2021

Signature/Date