

Product Number:	100457
Lot Number:	127514
Expiration:	2-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 24 Dec 2021

Signature/Date