

Product Number:	100457
Lot Number:	024112
Expiration:	28-Aug-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 05 Mar 2021

Signature/Date