

Product Number:	100453
Lot Number:	108831
Expiration:	29-Mar-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 15 Jun 2021
Signature/Date