

Product Number:	101359
Lot Number:	108123
Expiration:	22-Mar-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 26 Mar 2021
Signature/Date