

Product Number:	103158
Lot Number:	108008
Expiration:	21-Mar-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 24 Mar 2021
Signature/Date