

Product Number:	103158
Lot Number:	131411
Expiration:	10-Nov-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 08 Dec 2021
Signature/Date