

Product Number:	103158
Lot Number:	111221
Expiration:	22-Apr-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 05 May 2021
Signature/Date