

Product Number:	101362
Lot Number:	133915
Expiration:	5-Dec-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Sruvani Goel
Name

QC
Title

Sruvani Goel 18Jan2022
Signature/Date