

Product Number:	101956
Lot Number:	124613
Expiration:	3-Sep-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 15 Sep 2021

Signature/Date