

Product Number:	101359
Lot Number:	126809
Expiration:	25-Sep-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 13Oct2021
Signature/Date