

Product Number:	101955
Lot Number:	124910
Expiration:	6-Sep-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

Quality Control
Title

Shivani Goel 15 Sep 2021
Signature/Date