

Product Number:	101955
Lot Number:	204610
Expiration:	15-Feb-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

Quality Control
Title

Shivani Goel 24Feb2022
Signature/Date