

Product Number:	101955
Lot Number:	120410
Expiration:	23-Jul-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shwam Goel
Name

Quality Control
Title

Shwam Goel 23 Aug 2021
Signature/Date