

Product Number:	103158
Lot Number:	120030
Expiration:	19-Jul-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shwam Goel

Name

QC

Title

Shwam Goel 04 Aug 2021

Signature/Date