

Product Number:	100376
Lot Number:	120029
Expiration:	19-Jul-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 29 Jul 2021

Signature/Date