

Product Number:	100376
Lot Number:	112808
Expiration:	8-May-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 01Jun2021
Signature/Date