

Product Number:	100376
Lot Number:	104727
Expiration:	16-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 04 May 2021
Signature/Date