

Product Number:	100457
Lot Number:	109014
Expiration:	31-Mar-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 14 Apr 2021
Signature/Date