

Product Number:	100376
Lot Number:	133217
Expiration:	28-Nov-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval Shivani Goel QC Shivani Goel 18Jan2022
Name Title Signature/Date