

Product Number:	100376
Lot Number:	128906
Expiration:	16-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 01 Nov 2021

Signature/Date