

Product Number:	100457
Lot Number:	127518
Expiration:	2-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 15 Feb 2022
Signature/Date