

Product Number:	100457
Lot Number:	110524
Expiration:	15-Apr-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 17 Aug 2021
Signature/Date