

Product Number:	100453
Lot Number:	127411
Expiration:	1-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 12Oct2021
Signature/Date