

Product Number:	100458
Lot Number:	135012
Expiration:	16-Dec-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 15Feb2022
Signature/Date