

Product Number:	100457
Lot Number:	127516
Expiration:	2-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shwani Goel
Name

QC
Title

Shwani Goel 18 Jan 2022
Signature/Date