

Product Number:	101355
Lot Number:	127517
Expiration:	2-Oct-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 18Jan2022
Signature/Date