

Product Number:	101352
Lot Number:	122511
Expiration:	13-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 13 Oct 2021
Signature/Date