

Product Number:	100458
Lot Number:	105417
Expiration:	23-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 22 Mar 2021
Signature/Date