

Product Number:	100458
Lot Number:	122710
Expiration:	15-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 06 Oct 2021
Signature/Date