

Product Number:	101362
Lot Number:	934018
Expiration:	6-Dec-2020
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 06 Apr 2022
Signature/Date