

Product Number:	101362
Lot Number:	210305
Expiration:	13-Apr-2023
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 20 Apr 2022
Signature/Date