

Product Number:	102908
Lot Number:	205702
Expiration:	26-Feb-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 07 Mar 2022
Signature/Date