

Product Number:	102908
Lot Number:	127407
Expiration:	1-Oct-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.90-8.10

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 06Oct2021
Signature/Date