

Product Number:	101476
Lot Number:	134401
Expiration:	10-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95-9.05

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 15 Dec 2021
Signature/Date