

Product Number:	101475
Lot Number:	135005
Expiration:	16-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 06 Apr 2022
Signature/Date