

Product Number:	100460
Lot Number:	203013
Expiration:	30-Jan-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 02Feb2022
Signature/Date