

Product Number:	101358
Lot Number:	120910
Expiration:	27-Jul-2022
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 24 Dec 2021
Signature/Date