

Product Number:	101556
Lot Number:	113019
Expiration:	10-May-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 15 Dec 2021
Signature/Date