

Product Number:	103707
Lot Number:	120025
Expiration:	19-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 05 Aug 2021
Signature/Date