

Product Number:	103847
Lot Number:	211808
Expiration:	28-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval Shivani Goel QC Shivani Goel 03 May 2022
Name Title Signature/Date