

Product Number:	101353
Lot Number:	210412
Expiration:	14-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Shivani Goel QC manager
Richa Shrestha QC Supervisor Shwam Goel 03May2022
Name 03May2022 Title 03May2022 Signature/Date