

Product Number:	103360
Lot Number:	207610
Expiration:	17-Mar-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Shivani Goel
~~Richa Shrestha~~

Name

QC

Title

Shivani Goel 06 May 2022

Signature/Date