

Product Number:	103358
Lot Number:	105613
Expiration:	25-Feb-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Shivani Goel
Name

Quality Control *Manager*
Title

Shivani Goel 11Mar
Signature/Date 2021