

Product Number:	101476
Lot Number:	119011
Expiration:	9-Jul-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95-9.05

Review/Approval

Shivani Goel
~~Junpil Park~~

Manager, QC
~~Supervisor, QC~~

Shivani Goel 14 Jul 2021

Name

Title

Signature/Date