

Product Number:	102912
Lot Number:	210924
Expiration:	19-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Sonia Khillar

Name

QC

Title

Soni 12/1/29Aug202

Signature/Date