

Product Number:	101955
Lot Number:	229812
Expiration:	25-Oct-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Sonia Khilka

Name

QC Manager

Title

Sonia Khilka / 14 Nov 2022

Signature/Date