

Product Number:	101556
Lot Number:	223810
Expiration:	26-Aug-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 12/14/2023
Signature/Date