

Product Number:	101362
Lot Number:	316412
Control Number:	316412A
Expiration:	13-Jun-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khill 29 JUN 2023
Signature/Date