

Product Number:	101954
Lot Number:	312403
Control Number:	312403A
Expiration:	4-May-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sonia Killian
Name

QC Manager
Title

Sonia Killian 12MAY2023
Signature/Date