

Product Number:	101352
Lot Number:	302410
Expiration:	24-Jan-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Killian
Name

QC Manager
Title

Sonia Killian 1/27 JAN 2023
Signature/Date