

Product Number:	101362
Lot Number:	318008
Control Number:	318008A
Expiration:	29-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 Khillan 12 JUL 2023*

Signature/Date