

Product Number:	101362
Lot Number:	312908
Control Number:	312908A
Expiration:	9-May-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

[Signature] 23 MAY 2023
Signature/Date