

Product Number:	101476
Lot Number:	310202
Control Number:	310202A
Expiration:	12-Apr-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95 - 9.05

Review/Approval

Brian Ciccariello

Name

VP, QA+RA

Title

B-C 19 APR 2023

Signature/Date