

Product Number:	100376
Lot Number:	310014
Expiration:	10-Apr-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Brian Ciccarullo
Name

VP, QA/RA
Title

B - C *19 APR 2023*
Signature/Date