

Product Number:	101956
Lot Number:	316509
Control Number:	316509A
Expiration:	14-Jun-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 12 JUL 2023
Signature/Date