

Product Number:	101353
Lot Number:	313512
Control Number:	313512A
Expiration:	15-May-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 20 JUL 2023
Signature/Date