

Product Number:	101353
Lot Number:	319815
Control Number:	319815A
Expiration:	17-Jul-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 Sonia Khillan 21 Aug 2023
Title Signature/Date