

Product Number:	100453
Lot Number:	319116
Control Number:	319116A
Expiration:	10-Jul-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval Sonia Khillan QC Manager Sonia Khillan 20JUL2023
Name Title Signature/Date