

Product Number:	101357
Lot Number:	315711
Control Number:	315711A
Expiration:	6-Jun-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 29 JUN 2023
Name Title Signature/Date