

Product Number:	101357
Lot Number:	328314
Expiration:	10-Oct-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.18	6.18 - 6.22	PASS

Review/Approval

*Sonia Khillan*

Name

*Senior GC Analyst*

Title

*Sonia Khillan 19 Oct 2023*

Signature/Date