

Product Number:	101472
Lot Number:	329410
Expiration:	21-Oct-2024
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.35	7.35-7.45	PASS

Review/Approval

*Sonia Khiller*

Name

*Senior GC Analyst*

Title

*Sonia Khiller 12DEC2023*

Signature/Date