

Product Number:	103158
Lot Number:	334519
Expiration:	11-Dec-2024
Storage:	2-8 °C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.42	7.35 - 7.45	PASS

Review/Approval

Sonia Khillan
Name

Senior GC Analyst
Title

Sonia Khillan 21 DEC 2023
Signature/Date