

Product Number:	101472
Lot Number:	415510
Expiration:	3-Dec-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.39	7.35 - 7.45	PASS

Review/Approval

Robert Somogyi

QC Manager

 23 DEC 2024

Name

Title

Signature/Date