

Product Number:	101954
Lot Number:	416512
Expiration:	13-Dec-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.80	7.75 - 7.85	PASS

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 16.AUG.2024
Signature/Date