

Product Number:	103158
Lot Number:	413512
Expiration:	14-Nov-2025
Storage:	2-8°C

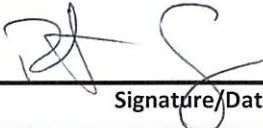
Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.37	7.35 - 7.45	PASS

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 26 JUN 2024
Signature/Date