

Product Number:	101476
Lot Number:	414115
Expiration:	20-May-2025
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	9.04	8.95 - 9.05	PASS

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 23MA/2024
Signature/Date