

Product Number:	101955
Lot Number:	414209
Expiration:	21-Nov-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.38	7.35 - 7.50	PASS

Review/Approval

Robert Somogyi

QC Manager

 13 JUN 2024

Name

Title

Signature/Date