

Product Number:	100457
Lot Number:	418314
Expiration:	1-Jan-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.35 - 7.45	PASS



Review/Approval

Robert Somogyi

QC Manager

Name

Title

  24 MAR 2025

Signature/Date