

Product Number:	101357
Lot Number:	505806
Control Lot Number:	505806A
Expiration:	27-Aug-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.19	6.18 - 6.22	PASS

Review/Approval

Robert Somogyi

QC Manager

 04 MAR 2025

Name

Title

Signature/Date