

Product Number:	100453
Lot Number:	505813
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

 27 MAR 2025

Name

Title

Signature/Date