

Product Number:	101357
Lot Number:	429524
Expiration:	21-Apr-2026
Storage:	2-8°C


**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	6.21	6.18 - 6.22	PASS

Review/Approval

Robert Somogyi

QC Manager

 22 JUN 2025

Name

Title

Signature/Date