

Product Number:	101357
Lot Number:	416609
Expiration:	14-Dec-2025
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

 28 JUN 2024

Name

Title

Signature/Date