

Product Number:	101352
Lot Number:	415103
Expiration:	30-Nov-2025
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

 05 JUN 2024

Name

Title

Signature/Date