

Product Number:	101474
Lot Number:	404514
Expiration:	14-Feb-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.35 - 7.45	PASS

RESULT

pH Reading	7.45
------------	------

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 16 FEB 2024

Signature/Date