

Product Number:	101362
Lot Number:	405807
Expiration:	27-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.75 - 7.85	PASS

RESULT

pH Reading	7.81
------------	------


Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 30 APR 2024

Signature/Date