

Product Number:	101474
Lot Number:	511201
Expiration:	22-Apr-2026
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.40
------------	------

Review/Approval

Robert Somogyi

QC Manager

Name

Title

Signature/Date

 23 APR 2025