

Product Number:	101359
Lot Number:	433808
Expiration:	3-Jun-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.30 - 7.50	PASS

RESULT

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

Review/Approval

Robert Somogyi
Name

Manager, Quality Control
Title


Signature/Date