

Product Number:	101359
Lot Number:	415512
Expiration:	3-Dec-2025
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.39
------------	------

Review/Approval

Robert Somogyi

QC Manager

Name

Title

 29 JAN 2025
Signature/Date