

Product Number:	100376
Lot Number:	426913
Expiration:	25-Mar-2026
Storage:	2-8°C

Laboratory Analysis

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

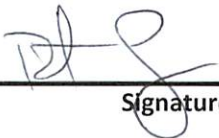
RESULT

pH Reading	7.37
------------	------

Review/Approval

Robert Somogyi

QC Manager

 09 OCT 2024

Name

Title

Signature/Date