

Product Number:	101359
Lot Number:	414210
Expiration:	21-Nov-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
Name

QC Manager
Title

 05 JUN 2024
Signature/Date