

Product Number:	101362
Lot Number:	405806
Expiration:	27-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.75 - 7.85	PASS

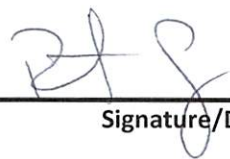
RESULT

pH Reading	7.81
------------	------

Review/Approval

ROBERT SOMOGYI  
Name

QC MANAGER  
Title

 16 APR 2024  
Signature/Date