

Product Number:	101474
Lot Number:	406707
Expiration:	7-Mar-2025
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.43
------------	------

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

 13 MAR 2024
Signature/Date