

Product Number:	103896
Lot Number:	411435
Expiration:	23-Oct-2025
Storage:	2 - 8°C

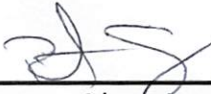
Laboratory Analysis

Characteristic	pH Target	Result	Disposition
pH	7.35 - 7.45	7.42	Pass

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 08 AUG 2024
Signature/Date