

Product Number:	104101
Lot Number:	406506
Expiration:	5-Sep-2025
Storage:	2 - 8°C

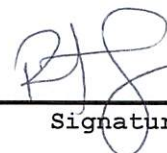
Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

 15 MAR 2024
Signature/Date