

Product Number:	104358
Lot Number:	324103
Expiration:	29-Oct-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

[Signature] 14 FEB 2024
Signature/Date