

Product Number:	104355
Lot Number:	326513
Expiration:	22-Nov-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

 12 SEP 2024
Signature/Date