

Product Number:	104101
Lot Number:	526709
Expiration:	24-Mar-2027
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title


25 Sep 2025
Signature/Date