

Product Number:	104101
Lot Number:	513906
Expiration:	19-Nov-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

<u>Herci Caswell</u>	<u>Director, QC</u>	<u>H. Caswell 03 June 2025</u>
Name	Title	Signature/Date