


Product Number:	102002
Lot Number:	405406
Expiration:	23-Aug-2025
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Robert Somogyi	QC Manager	 08 JAN 2025
Name	Title	Signature/Date