

Product Number:	101355
Lot Number:	415113
Expiration:	30-Nov-2025
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.35 - 7.45	PASS

Review/Approval

Binh Vu  
Name

QC Supervisor  
Title

 24 JUN 2024  
Signature/Date