

Product Number:	100457
Lot Number:	235405
Expiration:	20-Dec-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan  
Name

QC Manager  
Title

[Signature] / 03 JAN 2023  
Signature/Date