

Product Number:	102910
Lot Number:	320008
Expiration:	19-Jul-2024
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	Pass	7.95 - 8.05

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Kella 27 JUL 2023*  
Signature/Date