

Product Number:	100376
Lot Number:	120023
Expiration:	19-Jul-2022
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
pH Measurement	PASS

Review/Approval

*Shivani Goel*  
Name

QC  
Title

*Shivani Goel 23Jul2021*  
Signature/Date