

Product Number:	100376
Lot Number:	118214
Expiration:	1-Jul-2022
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 07Jul2021  
Signature/Date