

Product Number:	100458
Lot Number:	105414
Expiration:	23-Feb-2022
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 11 Mar 2021  
Signature/Date