

Product Number:	101355
Lot Number:	127507
Expiration:	2-Oct-2022
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Gopal  
Name

QC

Title

Shivani Gopal 30 Dec 2021

Signature/Date