

Product Number:	101355
Lot Number:	133218
Expiration:	28-Nov-2022
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel  
Name

QC

Title

Shivani Goel 02Feb2022  
Signature/Date