

Product Number:	101473
Lot Number:	121125
Expiration:	30-Jul-2022
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel  
Name

QC  
Title

Shwam Goel 26Jan 2022  
Signature/Date