

Product Number:	101476
Lot Number:	123306
Expiration:	21-Aug-2022
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	8.95-9.05

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 03 Sep 2021  
Signature/Date