

<b>Product Number:</b>	100460
<b>Lot Number:</b>	117224
<b>Expiration:</b>	21-Jun-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel  
Name

QC  
Title

Shwam Goel 13 Jul 2021  
Signature/Date