

<b>Product Number:</b>	101356
<b>Lot Number:</b>	304808
<b>Expiration:</b>	17-Feb-2024
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

<b>Review/Approval</b>	<i>Sonia Khilku</i> Name	<i>@C Manager</i> Title	<i>Sonia Khilku</i> 07 APR 2023 Signature/Date
------------------------	-----------------------------	----------------------------	---