

Product Number:	101353
Lot Number:	308307
Expiration:	24-Mar-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

<i>Sonia Khilkin</i>	<i>QC Manager</i>	<i>Sonia Khilkin</i> 07 APR 2023
Name	Title	Signature/Date