

Product Number:	101955
Lot Number:	327710
Expiration:	4-Oct-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.39	7.35 - 7.45	PASS

Review/Approval

<i>Sonia Khilkin</i>	<i>QC Manager</i>	<i>Sonia Khilkin 11 OCT 2023</i>
Name	Title	Signature/Date