

Product Number:	102910
Lot Number:	235005
Expiration:	16-Dec-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	Pass	7.95 - 8.05

Review/Approval	<u>Sonia Khilkin</u>	<u>QC Manager</u>	<u>Sonia Khilkin / 03 JAN 2023</u>
	Name	Title	Signature/Date