

Product Number:	101359
Lot Number:	234016
Expiration:	6-Dec-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval	<u>Sonia Khillan</u> Name	<u>QC Manager</u> Title	<u>Sonia KCH / 13DEC2022</u> Signature/Date
------------------------	------------------------------	----------------------------	------------------------------------------------