

Product Number:	101476
Lot Number:	326310
Expiration:	20-Sep-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.95	8.95 - 9.05	PASS

Review/Approval

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