

Product Number:	102914
Lot Number:	322311
Expiration:	11-Aug-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.98	7.95 - 8.05	PASS

Review/Approval

<i>Sonia Khillan</i>	<i>QC Manager</i>	<i>Sonia Khillan</i> 29 AUG 2023
Name	Title	Signature/Date