

Product Number:	101474
Lot Number:	322804
Expiration:	16-Aug-2024
Storage:	2 - 8 °C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.42	7.35 - 7.45	PASS

Review/Approval

Sonia Khillan QC Manager Sonia Kalla 01SEP2023

Name Title Signature/Date