

Product Number:	101474
Lot Number:	328202
Expiration:	9-Oct-2024
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia 12/11/2023
Signature/Date