

Product Number:	100376
Lot Number:	319209
Control Number:	319209A
Expiration:	11-Jul-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.37	7.35 - 7.45	PASS

Review/Approval

Sonia Khillan QC Manager
Name Title

Sonia Khillan 01SEP2023
Signature/Date