



Product Number:	101473
Lot Number:	319918
Control Number:	319918A
Expiration:	18-Jul-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.39	7.35 - 7.45	PASS

Review/Approval


Name

Title

Signature/Date

27 SEP 2023