

Product Number:	101472
Lot Number:	501525
Control Lot Number:	501525A
Expiration:	15-Jul-2026
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Robert Somogyi
Name

QC Manager
Title


Signature/Date