

Product Number:	101472
Lot Number:	501524
Control Lot Number:	501524A
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	QC Manager		29 JAN 2025
	Name	Title	Signature/Date	