

Product Number:	101356
Lot Number:	418313
Expiration:	1-Jan-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOSY
Name

QC MANAGER
Title

 14 NOV 2024
Signature/Date