

Product Number:	101909
Lot Number:	411503
Expiration:	24-Oct-2025
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] 06 MAY 2024
Signature/Date