

Product Number:	101473
Lot Number:	401622
Expiration:	16-Jul-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.38	7.35 - 7.45	PASS

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 24 SEP 2024
Signature/Date