

Product Number:	102914
Lot Number:	412407
Expiration:	3-May-2025
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.04	7.95 - 8.05	PASS

Review/Approval

ROBERT SOMOGYI

QC MANAGER

 15 MAY 2024

Name

Title

Signature/Date