

Product Number:	102912
Lot Number:	412405
Expiration:	3-May-2025
Storage:	2-8°C


Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.96	7.95 - 8.05	PASS

Review/Approval

ROBERT SOMOGYI

QC MANAGER

 15 MAY 2024

Name

Title

Signature/Date