

Product Number:	102908
Lot Number:	417606
Expiration:	24-Jun-2025
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.99	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi  
Name

QC Manager  
Title

 27 JUN 2024  
Signature/Date