

Product Number:	103847
Lot Number:	322712
Expiration:	15-Aug-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.42	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI

QC MANAGER

 20DEC2023

Name

Title

Signature/Date