

Product Number:	103358
Lot Number:	407910
Expiration:	19-Mar-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.02	7.95 - 8.05	PASS

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

 21 MAR 2024
Signature/Date