

Product Number:	103358
Lot Number:	419002
Expiration:	8-Jul-2025
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.00	7.95 - 8.05	PASS

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title


Signature/Date

23 JUL 2024