

Product Number:	103358
Lot Number:	431101
Expiration:	6-Nov-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.01	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 19 NOV 2024
Signature/Date