

Product Number:	103358
Lot Number:	505209
Expiration:	21-Feb-2026
Storage:	2 - 8°C

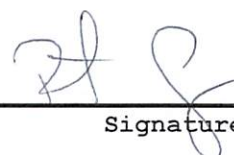
Laboratory Analysis

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

Review/Approval

Robert Somogyi

QC Manager

 25 FEB 2015

Name

Title

Signature/Date