

Product Number:	102908
Lot Number:	506415
Expiration:	5-Mar-2026
Storage:	2 - 8°C

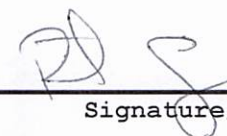
Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.96	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 10 MAR 2015
Signature/Date