

Product Number:	101956
Lot Number:	507617
Expiration:	17-Sep-2026
Storage:	2-8°C

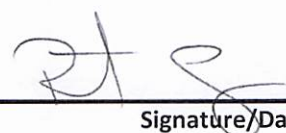
Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.35	7.35 - 7.45	PASS

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 21 MAR 2025
Signature/Date