

Product Number:	101476
Lot Number:	326309
Expiration:	20-Sep-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.95	8.95 - 9.05	PASS

Review/Approval

Pratish Shetty

Name

QC
SUPERVISOR

Title

Pratish Shetty

Signature/Date

27 SEP 2023