

Product Number:	101472
Lot Number:	324510
Expiration:	2-Sep-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.35 - 7.45	PASS

Review/Approval

Pratish S. Nit
Name

QC
Supervisor
Title

Pratish Nit
Signature/Date
27 SEP 2023