

Product Number:	101362
Lot Number:	333911
Expiration:	5-Dec-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.75 - 7.85	PASS

RESULT

pH Reading	7.85
------------	------

Review/Approval

Pratik Surti
Name

QC
SUPERVISOR
Title

Pratik Surti

08 DEC 2023

Signature/Date