

Product Number:	101474
Lot Number:	334013
Expiration:	6-Dec-2024
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.45
------------	------

Review/Approval

Pratish Jetti *QC* *Pratish Jetti* *08 DEC 2023*
Name Title Signature/Date