

Product Number:	101355
Lot Number:	304012
Expiration:	9-Feb-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia 12/1/24 FEB 2023
Signature/Date