

Product Number:	101473
Lot Number:	306119
Expiration:	2-Mar-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khilka QC Manager *[Signature]* 04APR2023
 Name Title Signature/Date