

Product Number:	101353
Lot Number:	234015
Expiration:	6-Dec-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Sonia Khilkin

Name

QC Manager

Title

Sonia Khilkin 08 DEC 2022

Signature/Date