

Product Number:	100453
Lot Number:	235403
Expiration:	20-Dec-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	6.18 - 6.22

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 103 JAN 2023

Signature/Date