

Product Number:	100457
Lot Number:	235406
Expiration:	20-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 10 FEB 2023
Signature/Date