

Product Number:	103360
Lot Number:	235009
Expiration:	16-Dec-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Sonia Khilka
Name

QC manager
Title

Sonia Khilka / 03JAN2023
Signature/Date