

Product Number:	100460
Lot Number:	405509
Expiration:	24-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

<u>ROBERT SOMOGYI</u>	<u>QC MANAGER</u>	<u>[Signature]</u> <u>04 MAR 2024</u>
Name	Title	Signature/Date