

Product Number:	103360
Lot Number:	412406
Expiration:	3-May-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.96	7.95 - 8.05	PASS

Review/Approval

<i>ROBERT SOMOGYI</i>	<i>QC MANAGER</i>	<i>[Signature]</i> 15 MAY 2024
Name	Title	Signature/Date