

Product Number:	101357
Lot Number:	404417
Expiration:	13-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.21	6.18 - 6.22	PASS

Review/Approval

Heidi Caswell

Director, QC

Heidi Caswell 14 Mar 2024

Name

Title

Signature/Date