

Product Number:	101357
Lot Number:	429523
Expiration:	21-Apr-2026
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 HCaswell 02Dec2024
Name Title Signature/Date