

Product Number:	101473
Lot Number:	426813
Expiration:	24-Mar-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.37	7.35 - 7.45	PASS

Review/Approval

Heidi Caswell
Name

Director, QC
Title

H. Caswell 26 Nov 2024
Signature/Date