

Product Number:	101352
Lot Number:	403115
Expiration:	31-Jul-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.44	7.35 - 7.45	PASS

Review/Approval

Heidi Caswell
Name

Director, QC
Title

Heidi Caswell 15 Feb 2024
Signature/Date