

Product Number:	101472
Lot Number:	514708
Expiration:	27-Nov-2026
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title

Kathleen Cornell
15 Aug 2025
Signature/Date