


Product Number:	101355
Lot Number:	428114
Expiration:	7-Apr-2026
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.40	7.35 - 7.45	PASS

Review/Approval      Robert Somogyi      QC Manager            05 DEC 2024

---

Name      Title      Signature/Date