

Product Number:	104072
Lot Number:	331019
Expiration:	6-May-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Heidi Caswell
Name

Director, IAC
Title

Heidi Caswell 14 Mar 2024
Signature/Date