

Product Number:	102002
Lot Number:	317904
Control Lot Number:	317904A
Expiration:	28-Jun-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

ROBERT SOMOGYI	QC MANAGER	 31 JAN 2024
Name	Title	Signature/Date