

Product Number:	100376
Lot Number:	108322
Expiration:	24-Mar-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Ruby Burke
Name

QC
Title

Ruby Burke 3/11/2021
Signature/Date