

Product Number:	100376
Lot Number:	104728
Expiration:	16-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Ruby Burke

Name

QC, Sr. Director

Title

Ruby Burke 05 Mar 2021

Signature/Date