

Product Number:	101556
Lot Number:	017801
Expiration:	26-Jun-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Patrizia Stadler

Associate Director, Quality
Control

P. Stadler 06 Jul 2020

Name

Title

Signature/Date