

Product Number:	101359
Lot Number:	002217
Expiration:	22-Jan-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Patrizia Stodler

Name

QC, Associate Director

Title

P. Stodler

30 Jun 2020

Signature/Date