

Product Number:	101359
Lot Number:	923111
Expiration:	19-Aug-2020
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Patrizia Stadler

Assoc. Director, QC

Name

Title

p. Stadler

19 Sep 2019

Signature/Date