

Product Number:	101357
Lot Number:	916503
Expiration:	14-Jun-2020
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler

19 Sep 2019

Signature/Date