



**G-Biosciences** ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** MOPS SDS Running Buffer [20X]

**Catalog No:** 786-927

**Release Date:** 11/5/19

**Size:** 1 gallon

**Retest Date:** 11/5/22

**Lot No:** 190511

**Storage:** Room temp.

### Specifications

**Appearance:** Clear, colorless to light yellow liquid

Test	Specification	Result
pH	7.60-8.51	7.95

This document has been produced electronically and is valid without signature

**Date of Print:** November 18, 2019

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)