



**G-Biosciences** ♦

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

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** MOPS Running Buffer [10X]

**Catalog No:** R031

**Release Date:** 11/5/20

**Size:** 1L

**Retest Date:** 11/5/23

**Lot No:** 200511

**Storage:** 4°C

### Specifications

**CAS Number:** N/A

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

Test	Specification	Result
pH	6.90-7.10	7.01

This document has been produced electronically and is valid without signature

**Date of Print:** November 13, 2020

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

Rev 10/25/18 BP



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