



**G-Biosciences** ♦

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

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** DMSO, Sterile Filtered

**Catalog No:** 786-1323 & 786-1388

**Release Date:** 1/30/20

**Size:** 10ml or 5 X 10ml

**Retest Date:** 1/30/24

**Lot No:** 203001

**Storage:** Room Temp.

### Specifications

**CAS Number:** 67-68-5

**Appearance:** Clear, colorless liquid

Test	Specification	Result
UV Absorbance 350nm	$\leq 0.01$	< 0.01
UV Absorbance 300nm	$\leq 0.10$	0.06
UV Absorbance 280nm	$\leq 0.30$	0.18
UV Absorbance 270nm	$\leq 0.70$	0.37
Purity (GC)	$\geq 99.70\%$	99.97%
Water (by Karl Fischer)	$\leq 0.2\%$	< 0.1%
Residue on Evaporation	$\leq 0.002\%$	< 0.001%

This document has been produced electronically and is valid without signature

**Date of Print:** February 5, 2020

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

Rev 10/25/18 BP



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