



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com



Certificate of Analysis

Product: Dimethylsulfoxide; DMSO

Catalog #: BKC-17

Lot #: 210906

Molecular Formula: $(\text{CH}_3)_2\text{SO}$

Molecular Weight: 78.13

Grade: Ultra Pure $\geq 99.0\%$

CAS# 67-68-5

<u>QC/QA Tests</u>	<u>Specification</u>	<u>Result</u>
Appearance	Clear Colorless Liquid	Passed
Protein NMR Spectrum	Consistent with Structure	Passed
Purity by Gas Chromatography	$\geq 99.0\%$ (Lot rejection if purity <99%)	99.5%

This document has been produced electronically and is valid without signature

For Research Use Only

Date of Print: October 12, 2021

G-Biosciences / Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Fax: 314-991-1504

DMSO C of A/Com # 3
Rev 04.19.13/IA



think proteins! think G-Biosciences!

www.GBiosciences.com