



Certificate of Analysis / Data Sheet

Product: DMSO, Sterile Filtered

Catalog No: 786-1322

Release Date: 8/7/20

Size: 1ml

Retest Date: 8/7/24

Lot No: 200708

Storage: -20°C

Specifications

CAS Number: 67-68-5

Appearance: Clear, colorless liquid

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

This document has been produced electronically and is valid without signature

Date of Print: August 18, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

