



BIOSCIENCES®

G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Glycine

Catalog No: RC-054

Size: 500g

Lot No: 200801

Release Date: 1/8/20

Retest Date: 1/8/23

Storage: Room Temp.

Specifications

CAS Number: 56-40-6

Appearance: White crystalline powder

Test	Specification	Result
Purity, dry matter	99.50-101.50%	100.0%
pH (5% in water)	5.8-6.2	6.2
Loss on Drying	≤ 0.1%	0.01%
Heavy Metals (Pb)	≤ 0.001%	< 0.001%
Arsenic (As)	≤ 0.0001%	< 0.001%
Residue on Ignition	≤ 0.08%	0.01%
Chloride	≤ 0.006%	< 0.006%
Sulfate (SO ₄)	≤ 0.006%	< 0.006%
Iron (Fe)	≤ 0.001%	< 0.001%
Lead (Pb)	≤ 0.0005%	< 0.0005%

This document has been produced electronically and is valid without signature

Date of Print: February 14, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com