



Certificate of Analysis / Data Sheet

Product: Guanidine Thiocyanate (Biotechnology Grade)
Catalog No: RC-059 **Release Date:** 9/2/20
Size: 500g **Retest Date:** 9/2/24
Lot No: 200209 **Storage:** Room Temp.

Specifications

CAS Number: 593-84-0

Appearance: White crystalline powder. Material may clump during storage, the quality of the product is not affected.

| Test | Specification | Result |
|-------------------------------|--------------------|---------------|
| Abs. @ 280nm (70%, water) | ≤ 0.8 | 0.3 |
| Abs. @ 300nm (6M, water) | < 0.10 | < 0.10 |
| Abs. @ 410nm (6M, water) | < 0.1 | 0.0084 |
| Copper | < 0.0001% | 0.000012% |
| DNase | None detected/Fail | None detected |
| Identification (IR) | PASS/Fail | PASS |
| Lead | < 0.0001% | 0.00004% |
| Melting Point | 118 - 121 °C | 119 °C |
| Protease | None detected/Fail | None detected |
| Purity | ≥ 99.0% | 100.0% |
| RNase | None detected/Fail | None detected |
| Solubility (70%, water 50 °C) | PASS/Fail | PASS |

This document has been produced electronically and is valid without signature

Date of Print: September 3, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

