



## Certificate of Analysis / Data Sheet

**Product:** Ellman's Reagent

**Catalog No:** BC87

**Release Date:** 11/1/19

**Size:** 5g

**Retest Date:** 11/1/23

**Lot No:** 190111

**Storage:** Room temp

### Specifications

**CAS Number:** 69-78-3

**Appearance:** Pale yellow powder

Test	Specification	Result
Melting Point	~241°C	241°C
Assay (aqueous acid-base titration)	≥98.5% to ≤101.5%	99.6%
Assay (HPLC)	≥98.5%	99.8%
Identification (FTIR)	Conforms	Conforms

This document has been produced electronically and is valid without signature

**Date of Print:** November 18, 2019

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

Rev 10/25/18 BP

