



## Certificate of Analysis / Data Sheet

**Product:** Urea

**Catalog No:** RC-112

**Size:** 1kg

**Lot No:** 201501

**Release Date:** 1/15/20

**Retest Date:** 1/15/26

**Storage:** Room Temp.

### Specifications

**CAS Number:** 57-13-6

**Appearance:** White crystalline powder

Test	Specification	Result
Identification (A, B)	Pass Test	PASS
Assay (Titration, dried basis)	99.0-100.5%	99.9%
Alcohol-Insoluble Matter	≤ 0.04%	0.01%
Chloride (Cl)	≤ 0.007%	0.0005%
Endotoxin	≤ 0.003 Units/mg	< 0.003 Units/mg
Heavy Metals (as Pb)	≤ 0.002%	0.0001%
Residual Solvents USP 467	Meets requirements	PASS
Residue on Ignition	≤ 0.1%	0.004%
Sterility	Meets requirements	PASS
Sulfate (SO <sub>4</sub> )	≤ 0.010%	0.0007%
Total impurities	≤ 2.0%	< 2.0%
Urea Related Compound A	≤ 0.1%	< 0.1%
BSE/TSE Free	BSE/TSE Free	PASS

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

