



Certificate of Analysis / Data Sheet

Product: Urea

Catalog No: RC-112

Release Date: 8/13/20

Size: 1kg

Retest Date: 8/13/26

Lot No: 201308

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%
Appearance of 20% solution	Pass Test	PASS
Residual Solvents USP 467	Meets requirements	PASS
Residue on Ignition	≤ 0.1%	0.1%
Individual Unspecified Impurities	≤ 0.1%	< 0.1%
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: September 2, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

