



Certificate of Analysis / Data Sheet

Product: Urea (Ultra Pure Grade)**Catalog No:** RC-112**Release Date:** 10/13/2021**Size:** 1kg**Retest Date:** 3/28/2022**Lot No:** 212209**Storage:** Room temp.**Specifications****CAS Number:** 57-13-6**Appearance:** White Crystalline Powder

Test	Specification	Result
Assay	99.0 – 100.5%	100.0%
Identification (A, B)	Pass Test	Passed
Appearance of 20% Solution	Pass Test	Passed
Alcohol- Insoluble Matter	≤ 0.04%	< 0.01%
Residue on Ignition	≤ 0.1%	< 0.01%
Urea Related Compound A	≤ 0.1%	< 0.1%
Individual Unspecified Impurities	≤ 0.1%	< 0.1%
Total impurities	≤ 2.0%	< 2.0%
Residual Solvents USP 467	Meets requirements	Passed
BSE/TSE Free	BSE/TSE Free	BSE/TSE Free
Grade	Ultra Pure	Ultra Pure
Country of Origin	Report	India

This document has been produced electronically and is valid without signature

Date of Print: January 28, 2022**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP

