



Certificate of Analysis / Data Sheet

Product: Guanidine Hydrochloride

Catalog No: BC85

Release Date: 2/4/20

Size: 250g

Retest Date: 2/4/24

Lot No: 200402

Storage: Room Temp.

Specifications

CAS Number: 50-01-1

Appearance: White crystal powder

Test	Specification	Result
pH	4.5-6.0	PASS
Loss on Drying	≤ 0.5%	PASS
Insoluble material	≤ 0.01% (in Water)	PASS
Related Substances	Ammonium chloride: ≤0.001%	PASS
Ash	≤ 0.05%	PASS
Assay by titration	≥99.70%	PASS

This document has been produced electronically and is valid without signature

Date of Print: October 9, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

