



Certificate of Analysis / Data Sheet

Product: Tris Base

Catalog No: RC-106

Size: 1kg

Lot No: 211509

Release Date: 9/23/21

Retest Date: 9/15/23

Storage: Room temp.

Specifications

CAS Number: 77-86-1

Appearance: White crystalline powder

Test	Specification	Result
Assay	99.5-101.0%	99.98%
MP	168-172°C	168.5-171.3°C
Insolubles	≤0.005%	Confirm
Heavy metals	≤5ppm	Confirm
Water content	≤0.5%	0.13%
pH (1.0% in water 25°C)	10.0-11.5pH unit	10.6
UV absorbance/260nm	≤0.1Abs unit	0.08
Fe	≤1ppm	Confirm
Cl	≤2ppm	Confirm

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

