



Certificate of Analysis / Data Sheet

Product: Trichloroacetic acid (TCA)

Catalog No: RC-813

Release Date: 10/13/21

Size: 500g

Retest Date: 9/21/23

Lot No: 212109

Storage: RT

Specifications

CAS Number: 76-03-9

Appearance: White to Off-White Powder or Crystal or Chunk(s)

Test	Specification	Result
Solubility (Color)	Colorless to Faint Yellow	Very Faint Yellow
Solubility (Turbidity) 1 g/ml, H ₂ O	Clear to Slightly Hazy	Clear
Infrared Spectrum	Conforms to Structure	Conforms
Titration with NaOH	≥ 99.0 %	99.8 %

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

