



Certificate of Analysis / Data Sheet

Product: Tris HCl

Catalog No: RC-107

Size: 500g

Lot No: 212307

Release Date: 7/28/21

Retest Date: 7/23/22

Storage: Room Temp.

Specifications

CAS Number: 1185-53-1

Appearance: White crystal powder

Test	Specification	Result
Assay(by Titn,dried basis)	Min.99.0%	99.5%
Melting Point(°C)	150.0-153.0°C	152.3
Loss on drying(105°C,3 hours)	≤0.5%	0.29%
Residue on Ignition	≤0.1%	0.02%
Solubility(1mlo/L Soln in H2O)	Clear	OK
Heavy metals Pb	≤5ppm	OK
Iron	≤1PPM	OK
pH(1.0% in water 25 °C)	4.2-5.0	4.6
pKa(20°C)	8.0-8.4	8.1
UVAbsorbance(260nm,1mol/L,1cm)	≤0.06	0.02
UV bsorbance(280nm,1mol/L,1cm)	≤0.05	0.009

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

