



Certificate of Analysis / Data Sheet

Product: EDTA, Disodium salt

Catalog No: 786-050

Release Date: 2/6/20

Size: 100g

Retest Date: 2/6/24

Lot No: 200602

Storage: Room Temp.

Specifications

CAS Number: 6381-92-6

Appearance: White crystalline powder

Test	Specification	Result
Assay (Dry Basis)	99.0-101.0%	99.75%
Identification (A,B,C)	Meets requirements	PASS
Solubility	Slightly soluble in water	PASS
pH (5%)	4.0-6.0	4.35
pH (1%)	4.3-4.7	4.48
Moisture (LOD)	8.7-11.4%	9.62%
Insoluble Matter	≤ 0.005%	0.003%
Nitroacetic Acid	≤ 0.1%	0.01%
Calcium	Conforms	PASS
Heavy Metals	≤ 50ppm	< 50ppm
Iron	≤ 0.01%	< 0.008%
Lead	≤ 10mg/kg	5mg/kg

This document has been produced electronically and is valid without signature

Date of Print: February 13, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

