



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** CTAB Extraction Buffer  
[2% CTAB, 100mM Tris, 1.4M NaCl, 20mM EDTA, pH 8.0]

**Catalog #:** 786-565

**Size:** 125ml

**Lot#:** 210408

**Appearance:** Clear liquid

### Tests:

Conductivity Test	PASSED
pH Test	PASSED
Specific gravity Test	PASSED

**Storage:** Store at 4°C.

*For Research Use Only*

This document has been produced electronically and is valid without signature

**Date of Print:** January 12, 2022

### **G-Biosciences / Genotech**

9800 Page Ave  
St. Louis, MO 63132, USA  
Fax: 314-991-1504

CTAB Extraction Buffer, 786-565 C of A/Com3  
Rev 12.21.10/IA



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)