



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

## Certificate of Analysis / Data Sheet

**Product:** TCEP-HCl; [Tris (2-carboxyethyl) phosphine, hydrochloride]  
**Catalog #:** 786-030  
**Size:** 1g  
**Lot #:** 182305  
**Grade:** Ultra Pure  
**Appearance:** White Solid  
**Molecular Formula:** C<sub>9</sub>H<sub>15</sub>O<sub>6</sub>P-HCl  
**Molecular Weight:** 286.7

<u>Tests</u>	<u>Specification</u>	<u>Result</u>
Solubility (1.1M)	Clear and colorless	Conforms
Loss on Drying	<1.0%	0.3%
Melting Point	177± 2°C	178°C
Purity (Assay)	≥ 98.0%	99.6%

**Storage:** Shipped at ambient temp, Store at -20°C, desiccated.

**Usage:** Used for the reduction of disulfide.

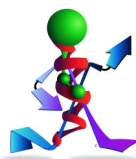
This document has been produced electronically and is valid without signature

**Date of Print:** November 2, 2018

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

9800 Page Avenue, St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504

TCEP-HCl C of A/Com # 3  
Rev 07.19.12/MM



think proteins! think G-Biosciences!

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