



Certificate of Analysis / Data Sheet

Product: TCEP-HCl; [Tris (2-carboxyethyl) phosphine, hydrochloride]
Catalog #: 786-030
Size: 1g
Lot #: 162904
Grade: Ultra Pure
Appearance: White Solid
Molecular Formula: C₉H₁₅O₆P-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: July 19, 2016

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

