



409PR-01

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

A Geno Technology, Inc. (USA) brand name

# Protein-S-S-Reductant™

(Cat. # 786-25PR)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Protein-S-S-Reductant™ contains neutralized and stabilized solution of 0.5M TCEP, suitable for reduction of protein before running electrophoresis - non-denaturing and SDS-PAGE denaturing. Protein-S-S-Reductant™ is a water soluble, odorless, non-toxic, and stable TCEP - Tri- (2-carboxyethyl) phosphine reductant for protein reduction. As compared to DTT and β-mercaptoethanol, the TCEP-reductant is more stable, more effective, and able to work over a wide range of pH, including lower acidic pH's. Reduces completely even the most stable disulfide bonds in less than 5 minutes at room temperature.

## ITEM(S) SUPPLIED (Cat# 786-25PR)

Description	Size
Protein-S-S-Reductant	4 x 1.0ml

## STORAGE CONDITIONS

It is shipped at ambient temperature. Store at -20°C, upon arrival. When stored and used properly, it is good for 1 year.

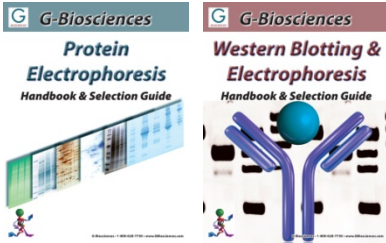
## PROTOCOL

### *Reduction of Disulfide bonds*

1. Add Protein-S-S-Reductant™ either in protein solution (20µl/ml) or 2X-electrophoresis sample loading buffer (40µl/ml).
2. After mixing with protein, vortex the tube and incubate the reaction for 5 minutes at room temperature.
3. At the end of incubation, the protein solution is ready for the next step. For SDS-PAGE, boil the protein sample for 5 minutes and load on gels.

## RELATED PRODUCTS

Download our Protein Electrophoresis and Western Blotting Handbooks.



<http://info.gbiosciences.com/complete-protein-electrophoresis-handbook/>

<http://info.gbiosciences.com/complete-western-blot-handbook--selection-guide/>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Last saved: 7/26/2012 CMH



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