



573PR-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

# SDS-PAGE Gel Fixing Solution

(Cat. #786-235, 786-236, 786-237)



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

## INTRODUCTION

The fixing of proteins in polyacrylamide gels prevents the diffusion of the proteins, thus keeping protein bands sharp and correctly resolved, and washes away running buffer components that may interfere with other downstream applications, such as staining. A major interfering agent is the SDS detergent used in SDS PAGE.

Protein fixation in polyacrylamide gels involves the denaturation and precipitation of the proteins into large insoluble aggregates within the gel matrix.

G-Biosciences SDS-PAGE Gel Fixing Solution is a relatively fast and convenient solution for fixing proteins in polyacrylamide gels and remove the SDS to prevent downstream interference.

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-235	SDS-PAGE Gel Fixing Solution	0.5L
786-236	SDS-PAGE Gel Fixing Solution	1L
786-237	SDS-PAGE Gel Fixing Solution	1gal

## STORAGE CONDITIONS

Shipped at ambient temperature. Upon receipt store at room temperature.

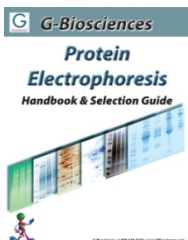
## PROTOCOL

1. Once protein electrophoresis is complete, transfer the protein gel to a suitable size container and cover the gel with the SDS-PAGE Gel Fixing Solution (~50ml)
2. Incubate for 1 hour at room temperature on an orbital shaker.
3. Discard the used SDS-PAGE Gel Fixing Solution and continue with downstream process.

**NOTE:** *Small or highly soluble proteins may not be sufficiently fixed by the above protocol, resulting in protein diffusion. We recommend prefixing with a 12% trichloroacetic acid solution for 1-3 hours.*

## RELATED PRODUCTS

Download our Protein Electrophoresis Handbook.



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

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

Last saved: 7/18/2013 CMH



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