



608PR-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

# PAGE-Optimizer™

For Preparing Samples for Gel Electrophoresis

(Cat. # 786-947, 786-948)



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

## INTRODUCTION

PAGE-Optimizer™ is a unique, spin column for the preparation of samples for SDS-PAGE. The columns contain a proprietary separation matrix that removes contaminants such as salts, detergents, cellular agents and other common buffering agents.

PAGE-Optimizer™ helps eliminate smeared lanes, distorted and/ or smiling bands and swollen lanes. PAGE-Optimizer™ has also been shown to reveal protein bands previously masked by the above interfering agents.

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-947	PAGE-Optimizer™	10 column
786-948	PAGE-Optimizer™	30 columns

## STORAGE CONDITIONS

The kit is shipped at ambient temperature. Store all the components at 4°C upon arrival.

## ADDITIONAL ITEMS REQUIRED

- Centrifuge Tubes
- Microfuge

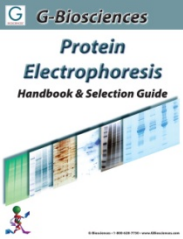
## PROTOCOLS

1. Before opening the column, centrifuge the PAGE-Optimizer™ for 30 seconds to collect the resin in the column.
2. Remove the cap and break open the bottom seal.
3. Place the PAGE-Optimizer™ in a 2ml collection tube and centrifuge at 1,000xg for 1 minute to remove the storage buffer. Discard the collection tube.
4. Position the PAGE-Optimizer™ in a clean 1.5ml tube and gently apply 5-10µl protein sample to the top of the resin bed.
5. Incubate at room temperature for 1 minute to allow the sample to enter the resin bed.
6. Centrifuge at 1,000xg for 5 minutes.
7. The sample collected in the tube is ready for SDS-PAGE analysis.

**NOTE:** PAGE-Optimizer™ columns are single use. For optimal results do not apply >10µl sample to a PAGE-Optimizer™ column. For processing larger sample volumes, use more than one column, each receiving a maximum of 10µl sample.

## RELATED PRODUCTS

Download our Protein Electrophoresis Handbook.



<http://info2.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: 3/9/2017 CMH



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