



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Gel Drying Solution

(Cat. # 786-684, 786-685, 786-686)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

G-Biosciences Gel Drying Solution effectively and evenly dries protein electrophoresis gels by the mechanism of passive evaporation and is suitable for permanent gel storage and densitometry etc. The Gel Drying Solution regulates the rate of drying and prevents the cracking of the gel. It is supplied in three convenient sizes; 500ml, 1L, and Gallon

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-684	Gel Drying Solution	500ml
786-685	Gel Drying Solution	1L
786-686	Gel Drying Solution	1gal

STORAGE CONDITIONS

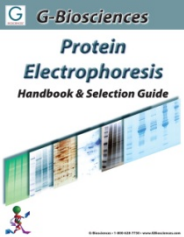
It is shipped at ambient temperature. Upon arrival, store it at room temperature. It is stable for one year when stored and used properly.

PROTOCOL

1. After staining and destaining the gel, wash three times with deionized water for 2-3 minutes each time on a rotary shaker.
2. Completely remove the water and add enough Gel Drying Solution to cover the gel
NOTE: For 8.5 x 7.5cm gel use approximately 35-40ml of Gel Drying Solution.
3. Equilibrate the gel in the Gel Drying Solution in a covered container, by shaking on a shaker for 15–20 minutes.
4. When equilibration process is complete, the gel is ready to be dried for permanent storage, densitometry etc.

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 or contact us.

Last saved: 2/14/2013 CMH



www.GBiosciences.com