



283PR

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

A Geno Technology, Inc. (USA) brand name

Geno ElectroPhore™

A Mini Horizontal Gel Electrophoresis System

(Cat. # MT-108)



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

ITEM(S) SUPPLIED:

1. Geno-ElectroPhore™ : A Mini Horizontal Gel Electrophoresis System (Cat # MT-108)

GENO-ELECTROPHORE™

Unpacking

Please verify that your MT-108 is complete with the following components.

- Electrophoresis tank with cover.
- Gel casting system (MT-G01)
- 1 Piece of 105 x 83 mm Tray (MT-T01)
- 2 Pieces of 50 x 83 mm Tray (MT-T02)
- 1 Piece of Analytical Comb 1.0 mm thick (MT-C01)
Two parts of 8 teeth for mini gel and one of 17 teeth for wide gel
- 1 Piece of Analytical Comb 1.0mm thick (MT-C02)
Two parts of 6 teeth for mini gel and one of 13 teeth for wide gel
- Power Cord (MT-P01)

Comb Dimension and Loading Volume Guide for MT-108

Cat. No.	Tooth Thickness (mm)	Tooth Width (mm)	Capacity per well ¹ (µl)
MT-C01	1.0	4.0	24
MT-C02	1.0	6.0	36

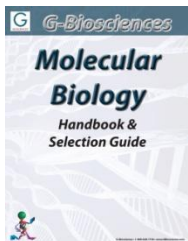
¹Capacity calculated for 6mm thick gel

Set-Up for Gel Casting

1. Put the Gel Tray in the casting system.
2. Select the analytical comb most suitable for your application.
3. Then pour 50-60ml of heated agarose gel into one of the 105 x 83mm trays or two of the 50 x 83mm trays.
4. Attach the comb to the comb holder.
5. Once the gel has formed, carefully pick up the comb and tray.

RELATED PRODUCTS

Download our Molecular Biology Handbook.



<http://info.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 6/16/2014 ADB



www.GBiosciences.com