



Geno-ElectroPhore™
A Mini Horizontal Gel Electrophoresis System
Cat # MT-108

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)

Table: Comb Dimension and Loading Volume Guide for MT-108

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

Capacity 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.

Note: For other related products, please visit our web site at www.GBiosciences.com
LU 11.04.08-SA/MM/IA

