

Introducing
Sequi-Gen GT Sequencing Cell

The Big Easy

GT Top Safety Cover
certified for IEC 1010-1
Safety Compliance

GT Lever Clamp
designed for easy gel casting

Upper buffer drainport
for easy buffer removal

GT Lever Clamp
attaches effortlessly and applies
uniform gel sandwich pressure

IPC (Integral Plate Chamber)
insures smile-free gels, everytime

GT Base
accommodates 4 different gel sizes

BIO-RAD
Sequi-Gen
Sequencing Cell



Casting gels with the Sequi-Gen GT systems is largely effortless.

The big, easy Sequi-Gen GT sequencing cell

The new Sequi-Gen GT cell is a big improvement over other electrophoresis systems. Its advanced design provides you the easiest way to manually cast and run nucleic acid sequencing gels. It includes many features that insure frustration-free usage, and smile-free results. In addition to IEC 1010-1 safety certification, the Sequi-Gen GT cell introduces a horizontal casting method that dramatically simplifies the process of gel casting. If you are looking for an easy gel casting method, this is big news.

Smile-free results deserve big grins

The Sequi-Gen GT systems give you smile-free results every time with our innovative Integral Plate Chamber (IPC). The IPC is a glass plate and plastic panel bonded together to form a thin (0.5 cm) chamber. This upper buffer chamber covers the entire area of the gel during electrophoresis. Heat generated during the run is quickly and evenly distributed across the gel by the buffer. Thus, temperature gradients causing sample mobility artifacts are eliminated. The IPC is easy to fill, easy to empty, and gives you results well worth grinning about.

This modular system is a big deal

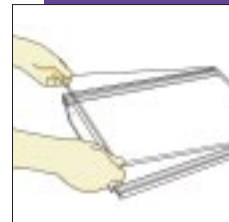
The Sequi-Gen GT base easily accommodates four different IPC sizes (21 x 40 cm, 21 x 50 cm, 38 x 30 cm, and 38 x 50 cm). The new 38 x 30 cm IPC is wide enough to load 100 samples and is designed for applications that require faster run times.

A huge choice of comb and spacer options

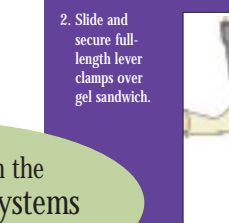
Many sharktooth and well-forming comb options are available for the Sequi-Gen GT cells. All combs and spacers are color-coded (based on thickness) and made from durable vinyl or plastic. Ask about our new multichannel pipet/microplate-compatible combs for high-throughput applications. They're sure to be immensely popular. For more information call 1-800-4BIORAD.

Sequi-Gen GT systems simplify a vast number of other applications

- Microsatellite analysis
- DNA fingerprinting
- Single-Strand Conformation Polymorphism (SSCP) studies
- Heteroduplex analysis
- DNA footprinting
- RNase protection assays
- S1 nuclease mapping
- Primer extension analysis
- DNA/protein binding studies (gel shift assays)
- Oligonucleotide analysis



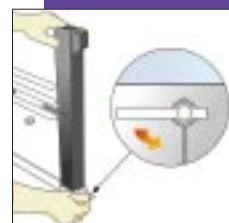
1. Assemble gel sandwich with IPC, outer glass plate, and spacers.



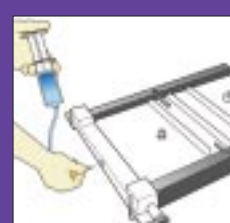
2. Slide and secure full-length lever clamps over gel sandwich.



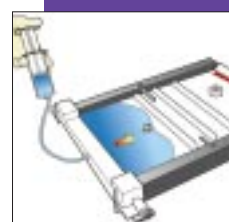
3. Insert gel sandwich assembly into precision caster base.



4. Lay assembly flat on lab bench and attach syringe tubing.



5. Inject gel solution between glass plates.



6. Position comb for gel polymerization.



Tape, grease, and binder clamps are eliminated. Gel casting does not get easier than with the new Sequi-Gen GT systems.

Ordering Information

Catalog Number	Product Description	Price	Catalog Number	Product Description	Price	Catalog Number	Product Description	Price
165-3860	Sequi-Gen GT System, 21 x 40 cm*	\$850.00	165-3802	Sequi-Gen GT/PowerPac 3000 System, 21 x 40 cm, 100/120 V [†]	2,800.00	165-3810	Sequi-Gen GT/PowerPac 3000 System, 38 x 30 cm, 100/120 V [†]	2,875.00
165-3861	Sequi-Gen GT System, 21 x 50 cm*	890.00	165-3805	Sequi-Gen GT/PowerPac 3000 System, 21 x 40 cm, 220/240 V [†]	Inquire	165-3811	Sequi-Gen GT/PowerPac 3000 System, 38 x 30 cm, 220/240 V [†]	Inquire
165-3862	Sequi-Gen GT System, 38 x 30 cm*	925.00	165-3803	Sequi-Gen GT/PowerPac 3000 System, 21 x 50 cm, 100/120 V [†]	2,840.00	165-3804	Sequi-Gen GT/PowerPac 3000 System, 38 x 50 cm, 100/120 V [†]	2,910.00
165-3863	Sequi-Gen GT System, 38 x 50 cm*	970.00	165-3806	Sequi-Gen GT/PowerPac 3000 System, 21 x 50 cm, 220/240 V [†]	Inquire	165-3807	Sequi-Gen GT/PowerPac 3000 System, 38 x 50 cm, 220/240 V [†]	Inquire

* Sequi-Gen GT systems include GT IPC assembly (IPC and bonded inner glass plate, outer glass plate, and clamp set), GT universal base, GT safety covers with cables, stabilizer bar, precision caster assembly (precision caster base, gasket, tubing, luer tapers, tubing, and syringes), 0.40 mm vinyl sharktooth comb and spacers, gel temperature indicator, leveling bubble, drain port connector, and instruction manual.

[†] Sequi-Gen GT/PowerPac 3000 systems include the appropriate Sequi-Gen GT system, PowerPac 3000 Power Supply, PowerPac temperature probe, and PowerPac instruction manual.

