

{ Just Add Water }



H Y D R O T E C H

A d v a n c e d G e l D r y i n g

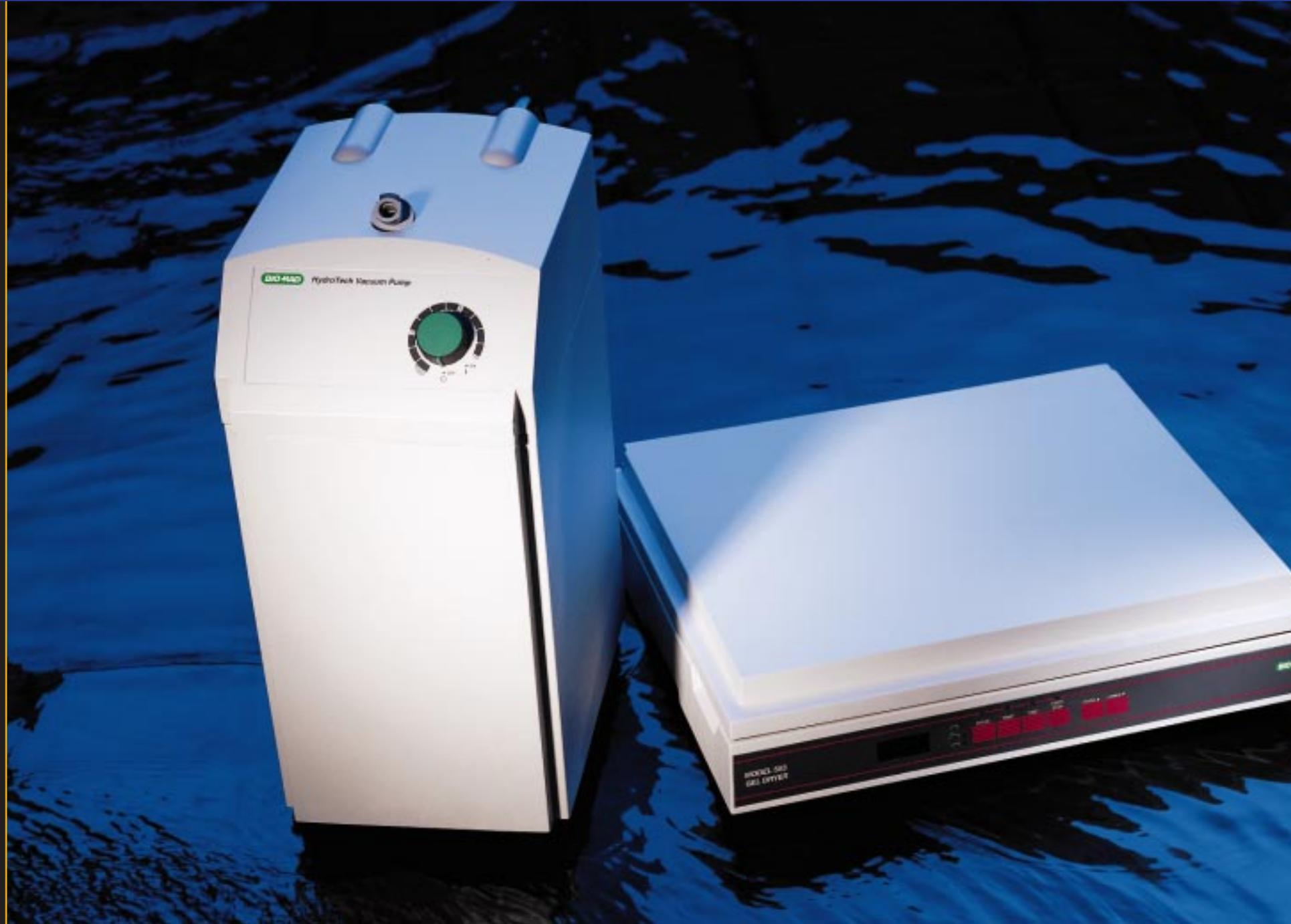
HYDROTECH

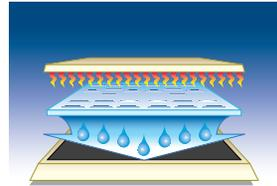
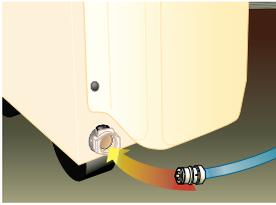
Gel Drying with cool, fresh water?

Introducing the new HydroTech Gel Drying System, a system that uses advanced, water-based vacuum technology to dry electrophoresis gels. It keeps the entire drying process refreshingly clean, pure and sensible.

The new HydroTech Vacuum Pump is packed with benefits:

- ◆ **NO oil changes, oil disposal or oil odor!**
The HydroTech is water-based. Say goodbye to oil pump maintenance, oil disposal costs and unpleasant oil odors.
- ◆ **NO liquid trap or vapor trap maintenance!**
All gel drying liquids and vapors are trapped in the water reservoir. Say goodbye to offensive chemical odors and the agony of trap maintenance.
- ◆ **NO expensive parts that continually need replacing!**
The HydroTech doesn't have expensive parts that wear out with normal use. Say goodbye to frequent and costly diaphragm replacements during the life of the pump.
- ◆ **Corrosion Resistant!**
The HydroTech won't corrode in the presence of common gel drying reagents. This may be the last gel drying pump you will ever need.
- ◆ **NO maintenance!**
{Just Add Water}. Emptying and refilling the water reservoir is the only maintenance you'll ever perform.





The Model 583 Gel Dryer

The HydroTech pump is designed to team up with the Model 583 Gel Dryer, the industry standard for efficiency and reliability.

◆ Precision Control

The Model 583 is microprocessor controlled which regulates temperature precisely and effortlessly.

◆ Exceptional Performance

The floating heating element heats gels from the top while the vacuum pulls liquid from the bottom. This unique top-down thermal flow produces uniform heating which prevents cracking and delivers unparalleled drying performance.

◆ Programmable

Three pre-programmed gel drying cycles to choose from: *Sequencing*, *Gradient*, *PAGE*.

Simply touch the keypad to select a pre-programmed drying cycle or program your own custom cycle.

◆ Easy Sealing

The quick seal track guarantees easy sealing of the transparent gasket. No more stretching or pulling of the gasket to seal the dryer. Closing the lid immediately seals the dryer.

Specifications

Model 583 Gel Dryer

Power requirements	100/120 volt models: 100-120 V, 50/60 Hz 220/240 volt models: 220-240 V, 50/60 Hz
Operating Time	5 minutes to 9 hrs 55 minutes
Operating cycles	3 pre-programmed drying cycles (sequencing, gradient, PAGE) and custom programming
Operating Temperature	50-90° C
Control System	Microprocessor controlled temperature and time
Heating Element	Floating (uniform heating)
Unique Thermal Flow	(heat applied from top)
Vacuum Requirements	Minimum vacuum 25" Hg (125 Torr) Minimum Capacity 20 liters/min

HydroTech Vacuum Pump

Ultimate Vacuum	>29" Hg (23.4 Torr) After 10 hrs >28 "Hg (48.8 Torr)
Pump Capacity	60 Hz: > 28 liters/min 50 Hz: > 21 liters/min
Fluid Capacity	MIN: 4 liters MAX: 5 liters
Function Mode	Continuous or Timed operation
Timer Control	0 - 3 hrs fully adjustable
Operating Conditions	4 - 35° C
Dimensions	40 x 25 x 51 cm (D x W x H)
Weight (dry)	14 kg (30 lb)
Power Requirements	100/120 V models: 100-120V, 50/60 Hz 220/240 V models: 220-240V, 50/60 Hz

Ordering Information

Catalog Number

Product Description

Gel Dryers

165-1745	Model 583 Gel Dryer , 100/120 V, includes power cord, porous gel support, transparent sealing gasket, filter paper backing, cellophane membrane backing, sequencing gel filter paper, and instruction manual
165-1746	Model 583 Gel Dryer , 220/240 V, includes power cord, porous gel support, transparent sealing gasket, filter paper backing, cellophane membrane backing, sequencing gel filter paper, and instruction manual

Gel Drying Systems

165-1789	HydroTech Gel Drying System , 100/120 V, includes Model 583 Gel Dryer, HydroTech Vacuum Pump, tubing and connectors
165-1790	HydroTech Gel Drying System , 220/240 V, includes Model 583 Gel Dryer, HydroTech Vacuum Pump, tubing and connectors

Vacuum Pumps

165-1781	HydroTech Vacuum Pump , 100/120 V, includes Vacuum Pump, power cord, quick disconnect fittings for 1/4 inch and 3/8 inch ID vacuum tubing, vacuum tubing, drain tubing, and instructions
165-1782	HydroTech Vacuum Pump , 220/240 V, includes Vacuum Pump, power cord, quick disconnect fittings for 1/4 inch and 3/8 inch ID vacuum tubing, vacuum tubing, drain tubing, and instructions
165-1788	HydroTech Vacuum Gauge , includes fittings to connect to the HydroTech vacuum line