

# NEW

## Double-Up Gel Dryer System



### Save Bench Space with the Double-Up Gel Dryer System

The new *Double-Up Gel Dryer System* combines two Model 583 Gel Dryers, a HydroTech™ vacuum pump, and the new Double-Up Gel Dryer Rack. For years, the Model 583 Gel Dryer has been the best selling slab dryer available. When teamed with the oil-less HydroTech vacuum pump, the Model 583 Gel Dryer forms a powerful combination for exceptional gel drying results.

The *Double-Up Gel Dryer Rack* accommodates two gel dryers up to 60 x 50 cm (W x D). The bottom shelf is on interlocking glides which allow full extension and easy access to the gel dryer's surface. The top shelf is stationary, and when a gel dryer is placed on it, provides stability to the rack. The Double-Up Gel Dryer Rack is plumbed for vacuum, made of sturdy sheet metal, and arrives assembled. The rack can be ordered separately, or as a system including two Model 583 Gel Dryers and one HydroTech vacuum pump.

**BIO-RAD**

## Specifications

Double-Up Gel Dryer Rack	
Overall dimensions	68 x 50 x 21 cm (W x D x H)
Bottom shelf dimensions	60 x 50 cm (W x D)
Accommodates dryers sizes	Up to 60 x 50 x 11 cm (W x D x H)
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 hours 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 above gel
Vacuum requirement	Minimum vacuum 25 inches Hg (125 Torr)
Minimum capacity	20 liters per minute
Safety certification (220/240V)	EN61010-1
Regulatory compliance (220/240V)	CE
HydroTech Vacuum Pump	
Ultimate vacuum	>29 inches Hg (23.4 Torr)
After 10 hours	>28 inches Hg (48.8 Torr)
Pump capacity	@ 60 Hz: >28 liters/min; @ 50 Hz: > 21 liters/min
Fluid capacity	Minimum: 4 liters; Maximum: 5 liters
Function mode	Continuous or timed operation
Timer control	0–3 hours fully adjustable
Operating conditions	4–35 °C
Dimensions	40 x 25 x 51 cm (D x W x H)
Weight (dry)	14 kg
Power requirements	115 V models : 100-120 V, 50/60 Hz; 230 V models: 220-240 V, 50/60 Hz
Safety certification (230 V)	EN61010-1
Regulatory compliance (230 V)	CE

## Ordering Information

Catalog number	Product Description
165-1797	Double-Up Gel Dryer System, 100/120 V, includes 2 Model 583 Gel Dryers, HydroTech vacuum pump, and Double-Up Rack
165-1798	Double-Up Gel Dryer System, 220/240 V
165-1796	Double-Up Gel Dryer Rack

**BIO-RAD**

**Bio-Rad  
Laboratories**

Life Science  
Group

Website [www.bio-rad.com](http://www.bio-rad.com) U.S. (800) 4BIORAD Australia 02-9914-2800 Austria (1)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771  
China (86-10) 2046622 Denmark 39 17 9947 Finland 90 804 2200 France (1) 43 90 46 90 Germany 089 31884-0 Hong Kong 7893300  
India 91-11-461-0103 Israel 03 951 4127 Italy 02-21609.1 Japan 03-5811-6270 The Netherlands 0313 18-540666 New Zealand 09-443 3099  
Singapore (65) 272-9877 Spain (91) 661 70 85 Sweden 46 (0) 8 627 50 00 Switzerland 01-809 55 55 United Kingdom 0800 181134