

The logo for Diagenode, featuring the word "diagenode" in a white, lowercase, sans-serif font. Above the letters "i", "e", and "o" are several small white dots of varying sizes, arranged in a slightly curved pattern. To the left of the text are several overlapping circles in shades of red and orange, some solid and some semi-transparent.

diagenode

Innovating Epigenetic Solutions

# DiaMag Rotator

**USER MANUAL**

VERSION 1

■ Cat. No. VL-100-0001

## 1. General specifications

Electrical requirements	
115V model 230V model	115V ± 10% - 50/60Hz, 0.03A 230V ± 10% - 50/60Hz, 0.015A
Speed 115V/230V model	24 rpm / 20 rpm
Motion	Rotating and rocking at 20°, 45°, 60°
Maximum tube capacity	36 x 1.5 ml or 2.0 ml micro tubes
Dimensions WxLxH	213x102x126 mm
Included accessories	Rotisserie for 36 x 1.5/2.0 ml micro tubes, set of two plastic pegs
Weight	0.8 kg
Operating conditions	4°C to 65°C, up to 85% RH, non-condensing

## 2. Installation of the Rotator

### Unpacking

Before starting the installation, carefully examine the rotator.

- Open the box and carefully remove the padding. Check that the rotator has not been visibly damaged during transportation. Keep the packing material until you are sure that the rotator works properly.
- Carefully lift the device out of the box.
- Be sure that the following information is included on the serial label located on the bottom of the unit: Serial Number, Product Number, Electrical rating.
- Check that the power cord plug has a pin configuration that matches a wall socket according to the local standard.

Should any kind of damage have occurred during transportation, contact the carrier immediately. The carrier is responsible for shipping damage.

### Powering the rotator

Place the rotator on horizontal, level and stable surface. Connect the plug end of the cord to a grounded wall socket. Avoid lines on which powerful electric motors, refrigerators and similar devices are connected. Use the on/off switch to control power to the rotator.



**Use caution when plugging into a grounded wall socket!**

**Do not touch plug with wet hands!**

**Do not unplug by pulling on the cord!**

## 3. Instructions for use

### Loading the rotisserie

The rotisserie supplied with the unit is designed to hold standard 1.5 ml or 2.0 ml microcentrifuge tubes. Install the rotisserie rod by screwing the two elements together through the hole in the center of the rotisserie plate. Insert the tubes into the tube holders. Mount the rotisserie in the apparatus in either the vertical position, using the rotisserie rod, or the horizontal position, using the bar keys on the plate. The side opposite to the driving mechanism is flexible to allow inserting the rotisserie without tools. Insert the bar key on the rotisserie rod or plate into the drive mechanism. The other end of the rotisserie is suspended in the bushing located in the opposite side of the unit. Be sure to distribute tubes evenly across the rotisserie carrier to achieve smooth rotating motion.

**To create tumbling motion** load the tubes and mount the rotisserie so the tubes are held in a vertical position. Then rest the unit on its end making sure that it is supported by the black side legs. The unit can be rested on either end. Setting the unit on the end with the power switch will create more vigorous tumbling motion than when it is rested on the opposite end.

**To create rocking motion** insert the plastic pegs supplied with the unit (stored in the holes at the lower part of the housing) into the holes located at the upper side of the unit. There are three sets of holes that will make the platform rock at 20°, 45° or 60°. When not in use store the pegs in the holes at the lower part of the housing.

**Mixing plates.** The rotisseries can be used for mixing in microplates. Remove the rotisserie rod if it is installed. Seal the microplates completely to avoid leakage. Make sure the microplate lines up with locking dimples. Fasten the plate (or plates) with rubber bands or small bungee cords (not supplied with the unit). Maximum of two deep well plates can be mixed at a time. When rocking, it is possible to shake just one plate. For all other types of motion, two plates should be placed on the rotisserie for proper balance.

## 4. Routine maintenance / cleaning

- No routine maintenance or lubrication is required.
- Keeping the rotator clean will help to maintain it in proper working order

The outside housing can be cleaned with a soft, damp cloth when needed.

Do not use any aggressive or abrasive cleaners (acetone, nitro, polish etc.) This may permanently damage the rotator finish.



**Before cleaning the rotator, always unplug the power cord from the wall socket!**

**Do not use the rotator near water sources. Be sure that water doesn't get on the device, especially during the cleaning procedures!**

- Do not expose the rotator to aggressive vapors (e.g. organic solvents).
- Hazardous samples should be mixed only in appropriate containment vessels.

NOTE: Misuse or unauthorized modifications will waive the warranty rights. If the equipment is used in a manner not specified by the manufacturer, the safety features of the rotator may be impaired.

**Do not mix flammable or explosive samples!**

## 5. Troubleshooting guide / service



Be sure to disconnect the rotator from the main power source immediately in case of any signs of malfunction and contact Diagenode.

## Ordering information

Description	Cat. No.
DiaMag Rotator	VL-100-0001
<b>DiaMag anti-mouse IgG-coated magnetic beads</b>	
DiaMag anti-mouse IgG-coated magnetic beads	kch-821-220
DiaMag anti-mouse IgG-coated magnetic beads	kch-821-660
DiaMag anti-mouse IgG-coated magnetic beads	kch-821-150
<b>DiaMag protein A-coated magnetic beads</b>	
DiaMag protein A-coated magnetic beads	kch-802-005
DiaMag protein A-coated magnetic beads	kch-802-220
DiaMag protein A-coated magnetic beads	kch-802-660
DiaMag protein A-coated magnetic beads	kch-802-150
<b>DiaMag protein G-coated magnetic beads</b>	
DiaMag protein G-coated magnetic beads	kch-818-005
DiaMag protein G-coated magnetic beads	kch-818-220
DiaMag protein G-coated magnetic beads	kch-818-660
DiaMag protein G-coated magnetic beads	kch-818-150
<b>DiaMag - magnetic racks</b>	
DiaMag02 - magnetic rack	kch-816-001
DiaMag1.5 - magnetic rack & disc stand	kch-716-015



# diagenode

Innovating Epigenetic Solutions

## Diagenode s.a. BELGIUM | EUROPE

Avenue de l'hôpital, 1  
Tour GIGA, 3rd Floor  
4000 Liège, Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

## Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101  
Denville, NJ 07834 USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

## | UK

Wellington House, East Road, Room 101  
CB1 1BH Cambridge, UK  
Tel: +44 1223 911396  
Fax: +44 1223 451100

For a complete listing of Diagenode's international distributors visit:  
<http://www.diagenode.com/pages/distributors.html>  
For rest of the world, please contact Diagenode s.a.