

diagenode

Innovating Epigenetic Solutions

# DiaMag Products for Epigenetic assays

Magnetic separation made easy and efficient



## Full Solution for your Epigenetic Research

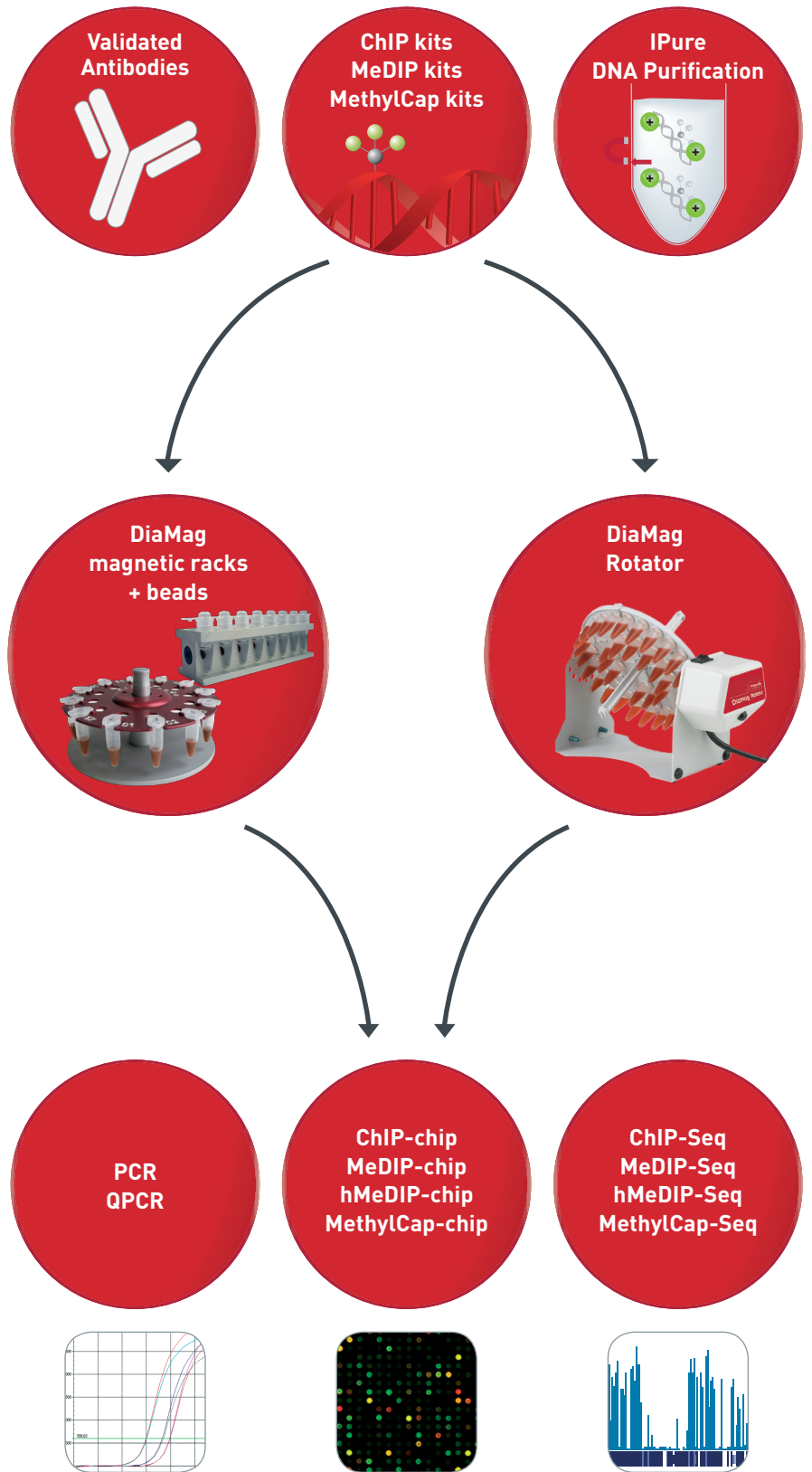
We offer a complete line of accessories and kits for ChIP and DNA methylation studies.

Diagenode's DiaMag products & technology allows you to get results quickly with easy handling and industry-leading reproducibility. Each DiaMag product fits seamlessly into your ChIP and MeDIP, hMeDIP, MethylCap and DNA Purification workflow and delivers the highest yields and quality of DNA recovery for any downstream application.

### The DiaMag Magnetic Bead Advantage

DiaMag Magnetic beads for immunoprecipitation provide numerous advantages over agarose beads as they allow for:

- Speed and versatility:** Small, uniform magnetic beads enable faster binding kinetics and can bind larger protein complexes unlike porous agarose beads.
- Higher specificity:** Magnetic beads have more specific binding and lower background due to less active binding potential.
- Greater sample yields and quicker workflow:** Magnetic beads do not require centrifugation reducing liquid handling, sample exposure to stress/destructive proteases, bead loss, and faster results.
- Cost efficiency:** Smaller amounts of magnetic beads are required as there is no centrifugation. In addition, scaling down bead amounts favors enrichment specificity.
- High throughput:** DiaMag magnetic beads are compatible with the IP-Star automation system.
- Application flexibility:** Magnetic beads allow for indirect ChIP with epitope tag-based approaches and bridge/secondary antibody binding capability.



## DiaMag Product Range - Features & Benefits

### DiaMag magnetic racks – fast, efficient isolation of magnetic beads in ChIP and MeDIP

- **High-throughput:** Processes 12-16 samples simultaneously
- **Standard:** Uses standard 1.5 and 0.2 ml tubes
- **High yields:** Prevents sample loss with efficient magnetic binding
- **Specific:** Eliminates sample contamination due to no tube handling
- **Sample integrity:** Maintains constant 4°C
- **Versatile:** Allows volume range of 25 ul-1 ml
- **Compatible** with all kits for ChIP, MeDIP, MBD, and DNA purification.

### DiaMag Magnetic Bead

- **Extensively validated** for ChIP, MeDIP, hMeDIP, MethylCap and IPure
- **Versatile** – Vast offering of antibody isotypes (see ordering table)
- **High binding** capacity of 3ug of antibody per 10 µl beads.

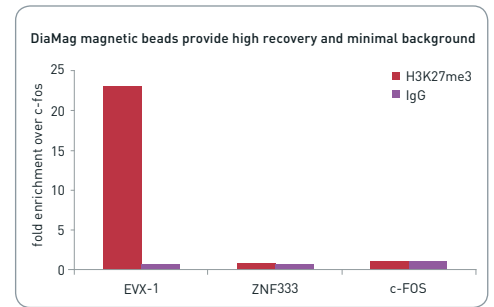
### DiaMag Rotator – seamless integration into your workflow

- **Compact:** Fits on your lab bench easily
- **Sample integrity:** Cold room compatible with gentle mixing
- **Standard:** Can process standard 1.5ml tubes and fits DiaMag 1.5 - magnetic rack tube holder.
- **High-throughput:** Holds up to 36 tubes of various sizes

### ChIP, MeDIP, MethylCap, MagBisulfite and IPure - magnetic bead-based kits

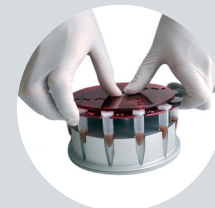
Obtain fast results with improved handling and reproducibility with our magnetic kits : Convenient kits with fewer steps, fewer buffers, and highly specific magnetic beads

- **HighCell# ChIP** – Recover large amounts of DNA from low abundance proteins without amplification and bias.
- **LowCell# ChIP** – Recover immunoprecipitated DNA in 24 hours with as few as 1,000 cells with this robust simplified kit
- **hMeDIP** – Obtain hydroxymethylation profiles in 24 hours with unique 5-hmC antibody and internal controls
- **MagMeDIP** - Obtain methylation profiles in 24 hours with unique 5-mC antibody and internal controls
- **MethylCap** – Obtain fractionation of methylated DNA by CpG density
- **MagBisulfite** – Obtain single base pair resolution of methylated DNA in less than 4 hours
- **IPure** – Delivers higher yields of pure, quality DNA for any downstream application after ChIP and MeDIP in comparison to standard column methods
- **Auto ChIP and DNA Methylation kits** – DiaMag magnetic beads are validated on Diagenode's IP-Star and allow for the complete automation of ChIP/MeDIP/hMeDIP/ and IPure.



ChIP results were obtained with the IP-Star® Automated System using the Auto ChIP kit containing DiaMag protein A-coated magnetic beads, which offer compatibility with rabbit polyclonal IgG antibodies. Ntera2 cells (100,000) were used per IP. Shown here is the enrichment of a positive genomic control (EVX-1) and two negative genomic controls (ZNF333 and c-FOS) using an H3K27me3 antibody.

### The DiaMag Racks and Rotator simplify your workflow without the need to work with individual tubes.



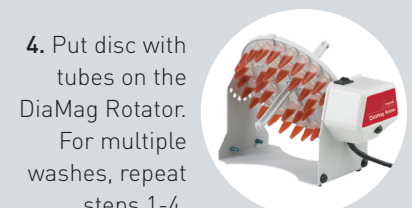
1. Place disc holder with tubes into rack. DiaMag beads adhere to the tube wall.



2. Remove the buffer with pipet without disrupting immobilized magnetic bead complex.



3. Take off disc with tubes and place onto disc stand. Resuspend sample with wash buffer.



4. Put disc with tubes on the DiaMag Rotator. For multiple washes, repeat steps 1-4.

Description	Cat. No.
<b>DiaMag magnetic racks and rotator</b>	
DiaMag02 - magnetic rack	kch-816-001
DiaMag1.5 magnetic rack & disc stand	kch-716-015
DiaMag1.5 - magnetic rack	kch-816-015
DiaMag1.5 - disc stand	kch-916-015
DiaMag Rotator	VL-100-0001
<b>DiaMag protein- or antibody-coated beads</b>	
DiaMag protein G-coated magnetic beads <i>Compatible with mouse IgG1, IgG2a, IgG2b and IgG3, rat IgG1, IgG2a, IgG2b and IgG3 rabbit and goat polyclonal Abs and human IgG1, IgG2, IgG3 and IgG4</i>	kch-818-220 (220 µl) kch-818-660 (660 µl) kch-818-150 (1.5 ml) kch-818-005 (5 ml)
DiaMag protein A-coated magnetic beads <i>Compatible with rabbit polyclonal Abs, mouse IgG2a, IgG2b and IgA, guinea pig IgG, dog IgG, pig IgG</i>	kch-802-220 (220 µl) kch-802-660 (660 µl) kch-802-150 (1.5 ml) kch-802-005 (5 ml)
DiaMag anti-mouse IgG-coated magnetic beads <i>Compatible with mouse IgG1, IgG2a, IgG2b</i>	kch-821-220 (220 µl) kch-821-660 (660 µl) kch-821-150 (1.5 ml)
<b>Magnetic bead-based kits</b>	
HighCell# ChIP kit	kch-mahigh-G16 (protein G, 16 rxns), kch-mahigh-A16 (protein A, 16 rxns)
LowCell# ChIP kit	kch-maglow-A16 (16 rxns), kch-maglow-A48 (48 rxns), kch-maglow-G16 (16 rxns), kch-maglow-G48(48 rxns)
hMeDIP kit* (monoclonal mouse antibody)	AF-110-0016
hMeDIP kit* (monoclonal rat antibody)	AF-104-0016
hMeDIP kit (polyclonal rabbit antibody)	AF-111-0016
MagMeDIP kit	mc-magme-A10 (10 rxns), mc-magme-048 (48 rxns)
MethylCap kit	AF-100-0048 (48 rxns)
IPure kit	AL-100-0100 (100 rxns)

\* Uses the only monoclonal antibody against 5-hydroxymethylcytosine (5-hmC) available in the market!

#### DIAGENODE HEADQUARTERS

#### 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

#### REGIONAL OFFICES

#### | 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.