

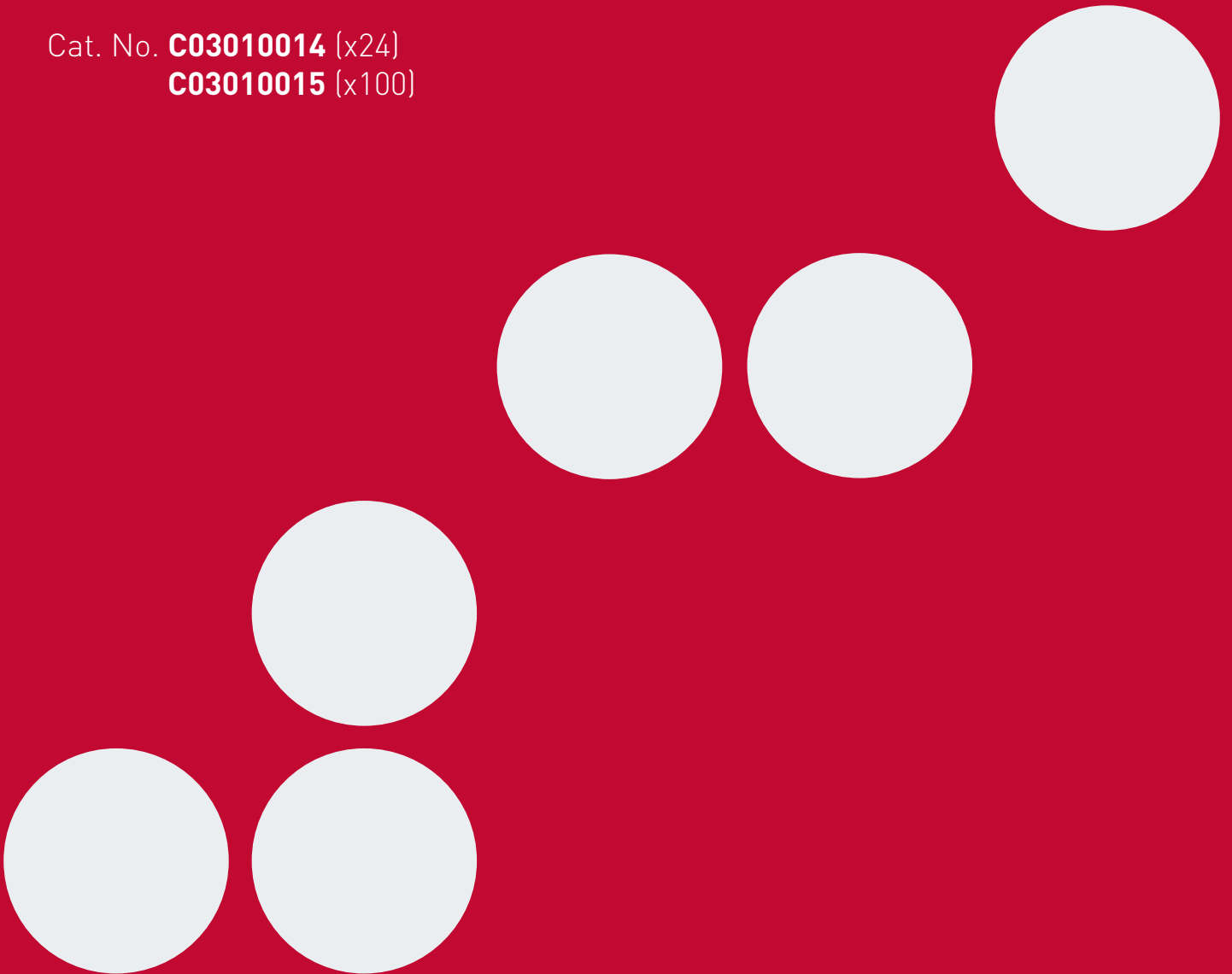


Innovating Epigenetic Solutions

# iPure kit v2

**Magnetic DNA Purification kit for epigenetic applications**

Cat. No. **C03010014** (x24)  
**C03010015** (x100)



# Ordering information

DIAGENODE HEADQUARTERS

## **Diagenode s.a. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing - 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  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

For a complete listing of Diagenode's international distributors visit:

<http://www.diagenode.com/company/distributors.php>

For rest of the world, please contact Diagenode sa.

Diagenode website: [www.diagenode.com](http://www.diagenode.com)

**FOR RESEARCH USE ONLY.**

**NOT INTENDED FOR ANY ANIMAL OR HUMAN THERAPEUTIC OR DIAGNOSTIC USE.**

© 2015 Diagenode SA. All rights reserved. No part of this publication may be reproduced, transmitted, transcribed, stored in retrieval systems, or translated into any language or computer language, in any form or by any means: electronic, mechanical, magnetic, optical, chemical, manual, or otherwise, without prior written permission from Diagenode SA (hereinafter, "Diagenode"). The information in this guide is subject to change without notice. Diagenode and/or its affiliates reserve the right to change products and services at any time to incorporate the latest technological developments. Although this guide has been prepared with every precaution to ensure accuracy, Diagenode and/or its affiliates assume no liability for any errors or omissions, nor for any damages resulting from the application or use of this information. Diagenode welcomes customer input on corrections and suggestions for improvement.

# Content

<b>Introduction</b> .....	<b>4</b>
<b>Method overview</b> .....	<b>5</b>
<b>Kit content</b> .....	<b>6</b>
<b>Time table</b> .....	<b>7</b>
<b>Protocol</b> .....	<b>8</b>
<b>Short protocol</b> .....	<b>8</b>
<b>Detailed protocol</b> .....	<b>10</b>
STEP 1a: Chromatin reverse crosslinking and elution (ChIP) .....	<b>10</b>
STEP 1b: DNA elution (MeDIP) .....	<b>11</b>
STEP 2: DNA binding (ChIP & MeDIP) .....	<b>12</b>
STEP 3: Washes (ChIP & MeDIP) .....	<b>12</b>
STEP 4: DNA elution (ChIP & MeDIP) .....	<b>13</b>
<b>Results</b> .....	<b>14</b>
<b>Diagenode's CHIP-seq workflow</b> .....	<b>15</b>

## Introduction

Diagenode's IPure kit is the only DNA purification kit using magnetic beads, that is specifically optimized for extracting DNA from ChIP and MeDIP (Chromatin IP and Methylated DNA IP).

It's a simple and straightforward protocol that delivers pure DNA ready for any downstream application (e.g. next generation sequencing). This approach guarantees a minimal loss of DNA and reaches significantly higher yields than a column purification (see results pag). Comparing to phenol-chloroform extraction, the IPure technology has the advantage of being nontoxic and much easier to be carried out on multiple samples. The use of the magnetic beads allows for a clear separation of DNA and increases therefore the reproducibility of your DNA purification.

**Two types of magnetic racks can be used:**



### **DiaMag1.5 - magnetic rack**

Cat. No. kch-716-015

- Holds 12x standard 1.5 ml tubes
- Working volume: 25-1000  $\mu$ l



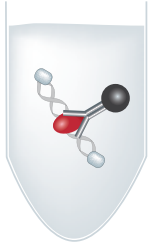
### **DiaMag02 - magnetic rack**

Cat. No. kch-816-001

- Holds 16x standard 0.2 ml tubes
- Working volume: 25-200  $\mu$ l

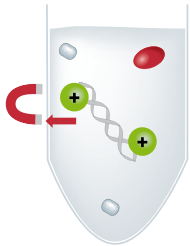
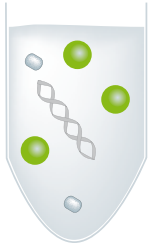
# Method overview

## IPure after ChIP



### STEP 1. Chromatin reverse cross-linking and elution

Chromatin is decrosslinked and eluted from beads (magnetic or agarose) which are discarded. **Magnetic beads** for purification are added.



### STEP 2. DNA binding

**Magnetic beads** acquire positive charge to bind the negatively charged phosphate backbone of DNA. DNA-bead complex is separated using a magnet.



### STEP 3. Washes

Proteins and remaining buffer are washed away.



### STEP 4. DNA Elution

DNA is eluted from magnetic beads, which are discarded. Purified DNA is ready for any downstream application:

- Next generation sequencing
- qPCR
- Amplification
- Microarray

## IPure after MeDIP



### STEP 1. DNA Elution

DNA is eluted from beads (magnetic or agarose). **Magnetic beads** for purification are added.



### STEP 2. DNA binding

**Magnetic Beads** acquire positive charge to bind the negatively charged phosphate backbone of DNA. DNA-bead complex is separated using a magnet.



### STEP 3. Washes

Remaining buffer is washed away.



### STEP 4. DNA Elution

DNA is eluted from magnetic beads, which are discarded. Purified DNA is ready for any downstream application:

- Next generation sequencing
- qPCR
- Amplification
- Microarray

## Kit content

The kit content is sufficient to perform 100 reactions.

IPure kit			
Description	Quantity (x24)	Quantity (x100)	Storage
200 µl tube strips (8 tubes/strip) + cap strips	4 pc	15 pc	Room temperature
Buffer A	2.4 ml	10 ml	4°C
Buffer B	96 µl	400 µl	4°C
Wash buffer 1 w/o isopropanol	1.5 ml	8 ml	4°C
Wash buffer 2 w/o isopropanol	1.5 ml	8 ml	4°C
Buffer C	1.2 ml	5 ml	4°C
IPure Beads v2	240 µl	1 ml	4°C
Carrier*	48 µl	200 µl	-20°C



\*This product is shipped at 4°C. Store it at -20°C upon arrival. This is an optimized buffer (NOT CARRIER DNA).

Related products			
Description	Old Cat. No.	New Cat. No.	Format
DiaMag02 - magnetic rack	kch-816-001	B04000001	1 pc
DiaMag1.5 - magnetic rack	kch-716-015	B04000003	1 pc
200 µl tube strips (8 tubes/strip) + cap strips	WA-002-0120	C30020002	120 pc
iDeal ChIP-seq kit for Histones x10	AB-001-0010	C01010050	10 rxns
iDeal ChIP-seq kit for Transcription Factors x10	/	C01010054	10 rxns
Auto iDeal ChIP-seq kit for Histones x24	/	C01010057	24 rxns
Auto iDeal ChIP-seq kit for Transcription Factors x24	/	C01010058	24 rxns
MagMeDIP kit x10	mc-magme-A10	C02010020	10 rxns
MagMeDIP kit x48	mc-magme-048	C02010021	48 rxns
Auto MeDIP kit x16	AF-Auto01-0016	C02010011	16 rxns
C02010012	AF-Auto01-0100	C02010012	100 rxns

### Required materials not provided

- DiaMag02 - magnetic rack (kch-816-001) / DiaMag1.5 - magnetic rack (kch-716-015)
- Microcentrifuge for 0.2 ml tubes or for 1.5 ml tubes with corresponding adaptor.
- 100% isopropanol
- Sodium Acetate
- 70% ethanol
- 100% ethanol
- DNase-free water

## Time table

IPure after ChIP (Chromatin Immunoprecipitation)		
STEP 1	Chromatin reverse crosslinking and elution	4 h (or overnight)
STEP 2	DNA binding	10 min
STEP 3	Washes	1 min
STEP 4	DNA elution	15 min
Total time		<b>4h30 min</b>

IPure after MeDIP (Methylated DNA Immunoprecipitation)		
STEP 1	DNA elution	30 min
STEP 2	DNA binding	10 min
STEP 3	Washes	1 min
STEP 4	DNA elution (option I / option II)*	15 min
Total time		<b>1h</b>

## Protocol

- » **Short**                      The 4 steps in 2 pages                      p.8
- » **Detailed**                      The 4 steps in 6 pages                      p.10

### Short protocol

After the last wash of immunoprecipitated material (ChIP or MeDIP), discard the last traces of wash buffer and use the pellet of beads (8-tube strip) for STEP 1a (ChIP) or STEP 1b (MeDIP).

#### STEP 1a: Chromatin reverse crosslinking and elution (ChIP)

1. Prepare the Elution buffer by mixing Buffer A and B as follows:

Elution buffer	1 rxn
Buffer A	115.4 µl
Buffer B	4.6 µl
<b>Total volume</b>	<b>120 µl</b>

2. Add 100 µl of Elution buffer to the bead pellets (tube strip).
3. Thaw your input sample (1.5 ml tube), and perform a pulse spin. Add 90 µl of Elution buffer and 10 µl of input sample to a new 200 µl tube (8 tube-strip).
4. Incubate samples and input for 4 hours (or overnight) at 65°C on a thermomixer, with continuous shaking.
5. Spin the 8-tube strip and place it into the DiaMag02 or DiaMag1.5 (magnetic rack). After 1 minute, transfer the supernatants to a new labelled 8-tube strip. Keep the samples on ice.

#### STEP 1b: DNA elution (MeDIP)

1. Prepare the Elution buffer by mixing Buffer A and B as follows:

Elution buffers	1 rxn
Buffer A	115.4 µl
Buffer B	4.6 µl
<b>Total volume</b>	<b>120 µl</b>

2. Add 50 µl of Elution buffer to the bead pellets (tube strip).
3. Thaw your input sample (1.5 ml tube), and perform a pulse spin. Add 92.5 µl of Elution buffer and 7.5 µl of input sample to a new 200 µl tube (8-tube strip).
4. Incubate samples and input DNA for 15 min at room temperature on a rotating wheel (40 rpm).
5. Spin the 8-tube strip and place it into the DiaMag02 or DiaMag1.5 (magnetic rack). After 1 minute, transfer the supernatants to a new labelled 8-tube strip.
6. Repeat the incubation of the bead pellets for 15 min at room temperature on a rotating wheel (40 rpm) in 50 µl elution buffer.
7. Spin the 8-tube strip. Place the 8-tube strips into the DiaMag02 or DiaMag1.5 containing the 50 µl IP samples, wait 1 minute and transfer the supernatants to the new labelled 8-tube strip to pool with the corresponding IP samples.



## STEP 2: DNA binding (CHIP & MeDIP)

1. Add **2 µl of carrier** to each IP and input sample. Vortex briefly and perform a short spin.
2. Add **100 µl of 100% isopropanol** to each IP and input sample. Vortex briefly and perform a short spin.



Following the addition of isopropanol the solution may become cloudy. This is not detrimental to your experiment and will not influence the quality or quantity of your purified DNA.

3. Resuspend the provided Magnetic beads and transfer 10 µl to each IP and input sample.
4. Incubate IP and input samples for 10 min at room temperature on a rotating wheel (40 rpm).

## STEP 3: Washes (CHIP & MeDIP)

1. Prepare the Wash buffer 1 containing 50% isopropanol for 100 reactions:

Wash buffer 1	24 rxns	100 rxns
Wash buffer 1 w/o isopropanol	1.5 ml	8 ml
Isopropanol (100%)	1.5 ml	8 ml
<b>Total volume</b>	<b>3 ml</b>	<b>16 ml</b>

2. Briefly spin the tubes, place into the DiaMag02 or DiaMag1.5, wait 1 minute and discard the buffer. Keep the captured beads and add per tube, 100 µl Wash buffer 1. Close the tubes, invert the 8-tube strip to resuspend the beads and incubate for 5 minutes at room temperature on a rotating wheel (40 rpm).
3. Prepare the Wash buffer 2 containing 50% isopropanol as follows:

Wash buffer 2	24 rxns	100 rxns
Wash buffer 2 w/o isopropanol	1.5 ml	8 ml
Isopropanol (100%)	1.5 ml	8 ml
<b>Total volume</b>	<b>3 ml</b>	<b>16 ml</b>

4. Briefly spin the tubes, place into the DiaMag02 or DiaMag1.5, wait 1 minute and discard the buffer. Keep the captured beads and add 100 µl Wash buffer 2 per tube. Close the tubes, invert the 8-tube strip to resuspend the beads and incubate for 5 minutes at room temperature on a rotating wheel (40 rpm).

## STEP 4: DNA elution (CHIP & MeDIP)

1. Briefly spin the tubes and place them into the DiaMag02 or DiaMag1.5, wait 1 minute and discard the buffer. Keep the captured beads and add per tube, 25 µl buffer C (alternatively, a higher volume can be used for the elution if necessary). Close the tubes, invert the 8-tube strip to resuspend the beads and incubate for 15 minutes at room temperature on a rotating wheel (40 rpm). Resuspend the pelleted beads using the pipet and make sure that you drop them on the bottom of the tube.
2. Spin the two 8-tube strips and place it into the DiaMag02 or DiaMag1.5, wait 1 minute and transfer the supernatants into a new labelled 1.5 ml tube. Keep the bead pellets on ice.
3. Place the DNA on ice and proceed to any desired downstream applications, or store it at -20°C or -80°C until further use.

## Detailed protocol

After the last wash of immunoprecipitated material (ChIP or MeDIP), discard the last traces of wash buffer and use the pellet of beads (8-tube strip) for STEP 1a (ChIP) or STEP 1b (MeDIP).

### STEP 1a. Chromatin reverse crosslinking and elution (ChIP)



1. Prepare the **Elution buffer** by mixing **Buffer A** and **B** as follows:

Elution buffer	1 rxn
Buffer A	115.4 $\mu$ l
Buffer B	4.6 $\mu$ l
<b>Total volume</b>	<b>120 <math>\mu</math>l</b>

- Place Buffer A at 25°C during 30 min before use.
- 100  $\mu$ l of Elution buffer are needed per IPure reaction (20  $\mu$ l excess).
- 1 IPure reaction corresponds to the purification 1 ChIP or 1 input sample.

#### Attention:

**Make sure when working with Buffer A, that there are no crystals left in solution. Otherwise heat up gently and mix until complete disappearance of such crystals.**

2. Add **100  $\mu$ l of Elution buffer** to the bead pellets (tube strip).
3. Thaw your input sample (1.5 ml tube), and perform a pulse spin. Add **90  $\mu$ l of Elution buffer** and **10  $\mu$ l of input sample** to a new 200  $\mu$ l tube (8 tube-strip).
  - Input sample corresponds to 10% of the IP sample.
4. Incubate samples and input for 4 hours (or overnight) at 65°C on a thermomixer, with continuous shaking.
5. Spin the 8-tube strip and place it into the **DiaMag02 or DiaMag1.5** (magnetic rack). After 1 minute, transfer the supernatants to a new labelled 8-tube strip. Keep the samples on ice.

## STEP 1b. DNA elution (MeDIP)



1. Prepare the **Elution buffer** by mixing **Buffer A** and **B** as follows:

Elution buffer	1 rxn
Buffer A	115.4 $\mu$ l
Buffer B	4.6 $\mu$ l
<b>Total volume</b>	<b>120 <math>\mu</math>l</b>

- Place Buffer A at 25°C during 30 min. before use.
- 100  $\mu$ l of Elution buffer are needed per IPure reaction (20  $\mu$ l excess).
- 1 IPure reaction corresponds to the purification 1 MeDIP or 1 input sample.

### Attention:

**Make sure when working with Buffer A, that there are no crystals left in solution. Otherwise heat up gently and mix until complete disappearance of such crystals.**

2. Add **50  $\mu$ l of Elution buffer** to the bead pellets (tube strip).
3. Thaw your input sample (1.5 ml tube), and perform a pulse spin. Add **92.5  $\mu$ l of Elution buffer** and **7.5  $\mu$ l of input sample** to a new 200  $\mu$ l tube (8-tube strip).
  - Input sample corresponds to 10% of the IP sample.
4. Incubate samples and input DNA for **15 min at room temperature** on a rotating wheel (40 rpm).
5. Spin the 8-tube strip and place it into the **DiaMag02 or DiaMag1.5** (magnetic rack). After 1 minute, transfer the supernatants to a new labelled 8-tube strip.
6. Repeat the incubation of the bead pellets for **15 min at room temperature** on a rotating wheel (40 rpm) in **50  $\mu$ l elution buffer**.
  - For input DNA samples: 1 elution in 100  $\mu$ l.
  - For MeDIP samples: 2 elutions in 50  $\mu$ l (total volume 100  $\mu$ l).
7. Spin the 8-tube strip. Place the 8-tube strips into the **DiaMag02 or DiaMag1.5** containing the 50  $\mu$ l IP samples, wait 1 minute and transfer the supernatants to the new labelled 8-tube strip to pool with the corresponding IP samples.
  - Elutions of IP and input samples are now completed in 100  $\mu$ l and are in the same 8-tube strip.

## STEP 2. DNA binding (ChIP & MeDIP)



1. Add **2 µl of carrier** to each IP and input sample. Vortex briefly and perform a short spin.
2. Add **100 µl of 100% isopropanol** to each IP and input sample. Vortex briefly and perform a short spin.



Following the addition of isopropanol the solution may become cloudy. This is not detrimental to your experiment and will not influence the quality or quantity of your purified DNA.

3. Resuspend the provided **Magnetic beads** and transfer 10 µl to each IP and input sample.
  - Keep Magnetic beads in liquid suspension during storage at 4°C and at all handling steps, as drying will result in reduced performance.
  - The final volume is now 212 µl per IPure reaction.
4. Incubate IP and input samples for 10 min at room temperature on a rotating wheel (40 rpm).

## STEP 3. Washes (ChIP & MeDIP)



1. Prepare the **Wash buffer 1** containing 50% isopropanol for 100 reactions:

Wash buffer 1	24 rxns	100 rxns
Wash buffer 1 w/o isopropanol	1.5 ml	8 ml
Isopropanol (100%)	1.5 ml	8 ml
<b>Total volume</b>	<b>3 ml</b>	<b>16 ml</b>

- Never leave the bottle open to avoid evaporation.
2. Briefly spin the tubes, place in the **DiaMag02 or DiaMag1.5**, wait 1 minute and discard the buffer. Keep the captured beads and add per tube, **100 µl Wash buffer 1**. Close the tubes, invert the 8-tube strip to resuspend the beads and incubate for 5 minutes at room temperature on a rotating wheel (40 rpm).
    - Do not disturb the captured beads attached to the tube wall.
    - Always briefly spin the tubes to bring down liquid caught in the lid prior to positioning into the Diagenode Magnetic Rack.
  3. Prepare the **Wash buffer 2** containing 50% isopropanol as follows:

Wash buffer 2	24 rxns	100 rxns
Wash buffer 2 w/o isopropanol	1.5 ml	8 ml
Isopropanol (100%)	1.5 ml	8 ml
<b>Total volume</b>	<b>3 ml</b>	<b>16 ml</b>

- Never leave the bottle open to avoid evaporation.
4. Wash the IP and input samples with the **Wash buffer 2** as follows. Briefly spin the tubes, place into the **DiaMag02 or DiaMag1.5**, wait 1 minute and discard the buffer. Keep the captured beads and add per tube, **100 µl Wash buffer 2**. Close the tubes, invert the 8-tube strip to resuspend the beads and incubate for 5 minutes at room temperature on a rotating wheel (40 rpm).
    - Do not disturb the captured beads attached to the tube wall.
    - Always briefly spin the tubes to bring down liquid caught in the lid prior to positioning in the **DiaMag02**.

## STEP 4. DNA elution (ChIP & MeDIP)



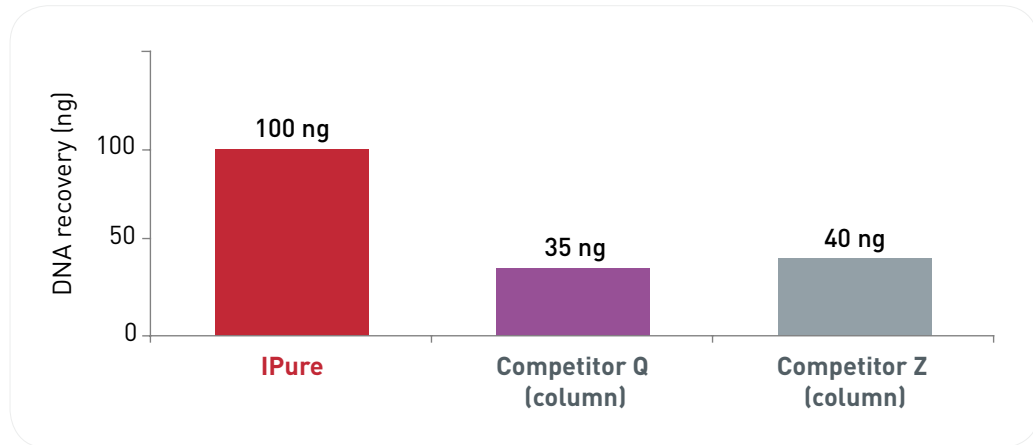
**Note:**

This elution buffer (Buffer C) is suitable for direct qPCR analysis, whole genome amplification, chip hybridization and next generation sequencing.

1. Briefly spin the tubes and place them into the **DiaMag02 or DiaMag1.5**, wait 1 minute and discard the buffer. Keep the captured beads and add per tube, **25 µl buffer C** (alternatively, a higher volume can be used for the elution if necessary). Close the tubes, invert the 8-tube strip to resuspend the beads and incubate for **15 minutes at room temperature** on a rotating wheel (40 rpm). Resuspend the pelleted beads using the pipet and make sure that you drop them on the bottom of the tube.
2. Spin the 8-tube strips and place it into the **DiaMag02 or DiaMag1.5**, wait 1 minute and transfer the supernatants into a new labelled 1.5 ml tube. Keep the bead pellets on ice.
3. Place the DNA on ice and proceed to any desired downstream applications, or store it at -20°C or -80°C until further use.

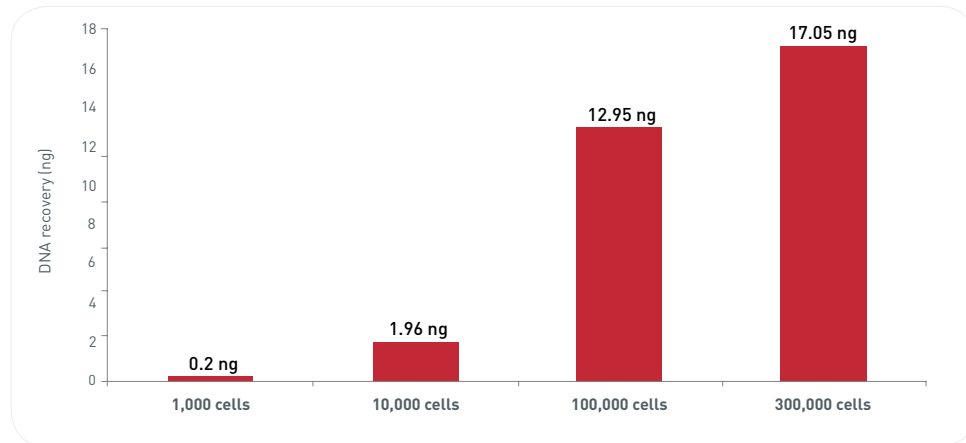
## Results

### Comparison of DNA recovery after purification with IPure technology and competitor kits



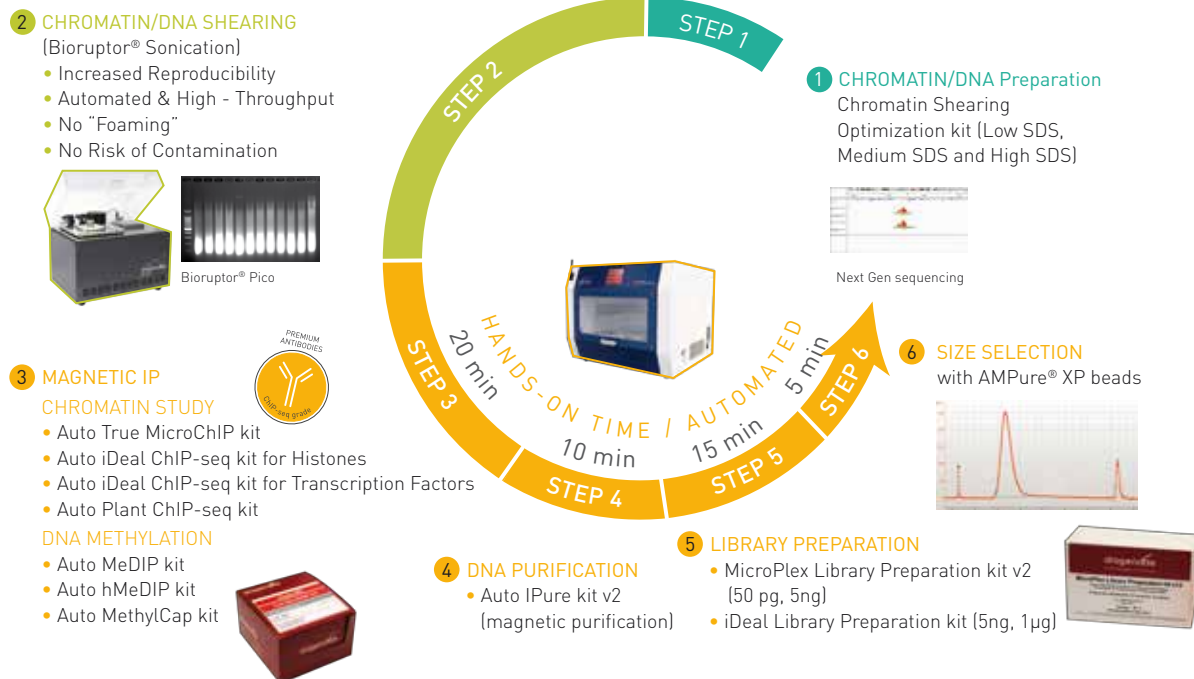
MeDIP assays were performed using the MagMeDIP kit (cat# mc-magme-048). The immunoprecipitated samples were purified with the IPure technology and two competitor kits (competitor Q and Z). The purified DNA was eluted in 50  $\mu$ l of water and quantified with a Nanodrop.

### DNA recovery after purification of ChIP samples using IPure technology



ChIP assays were performed using different amounts of U2OS cells, the LowCell# ChIP kit (cat# kch-maglow-016) and the H3K9me3 antibody (Cat. No. pAb-056-050; 2  $\mu$ g/IP). The purified DNA was eluted in 50  $\mu$ l of water and quantified with a Nanodrop.

## Bringing it all together: Diagenode's ChIP-seq workflow



**Figure 1. Diagenode provides a full suite of manual and automated solutions for ChIP experiments.**

For Step 1, we offer products to isolate nuclei and chromatin. Step 2 describes reproducible sample shearing with the Bioruptor® product line. In Step 3 and Step 4, the Diagenode IP-Star® Compact provides error-free, walk-away automation for all your immunoprecipitation and antibody capture needs.



Innovating Epigenetic Solutions

**DIAGENODE  
HEADQUARTERS**

**DIAGENODE S.A.  
BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing - 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  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**FOR A COMPLETE LISTING OF  
DIAGENODE'S INTERNATIONAL  
DISTRIBUTORS VISIT:**

<https://www.diagenode.com/>  
distributors For rest of the world,  
please contact Diagenode sa.  
[www.diagenode.com](http://www.diagenode.com)