



BIOARRAY™

methyated DNA IP Kit

Enrich for methylated DNA (100-fold) from sample for complete diverse downstream applications

The BIOARRAY™ methylated DNA IP kit features IP technology for the enrichment of 5-methylcytosine-containing DNA from any pool of fragmented genomic DNA for use in genome-wide methylation analysis. The kit includes a highly specific anti-5-methylcytosine monoclonal antibody for the “capture” and separation of methylated DNA from non-methylated DNA in only a few hours.

Methylation is a naturally occurring event in both prokaryotic and eukaryotic organisms. In prokaryotes DNA methylation provides a way to protect host DNA from digestion by restriction endonucleases that are designed to eliminate foreign DNA, and in higher eukaryotes DNA methylation functions in the regulation/control of gene expression.

Ordering Information

[Order Online »](#)

| | |
|----------------|----------|
| ENZ-45012-0010 | 10 tests |
|----------------|----------|

Manuals, SDS & CofA

[View Online »](#)

- Highly specific antibody is included to distinguish methylated DNA from pool
- DNA enrichment for large-scale DNA methylation analysis
- Easy to use with reproducible results

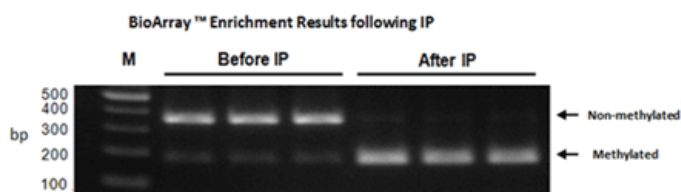


Figure demonstrates efficient enrichment of methylated vs. non methylated DNA by using the BioArray™DNA IP kit. Salmon sperm genomic DNA “spiked” with non methylated/methylated control DNA was processed with the kit. The eluted DNA was amplified by PCR using the supplied control primers. Digestion of the amplicons with Nco I produced two 175bp fragments for the methylated DNA control or one 350bp fragment for the non methylated DNA control. The results show an efficient enrichment of methylated versus non methylated DNA by the kit.

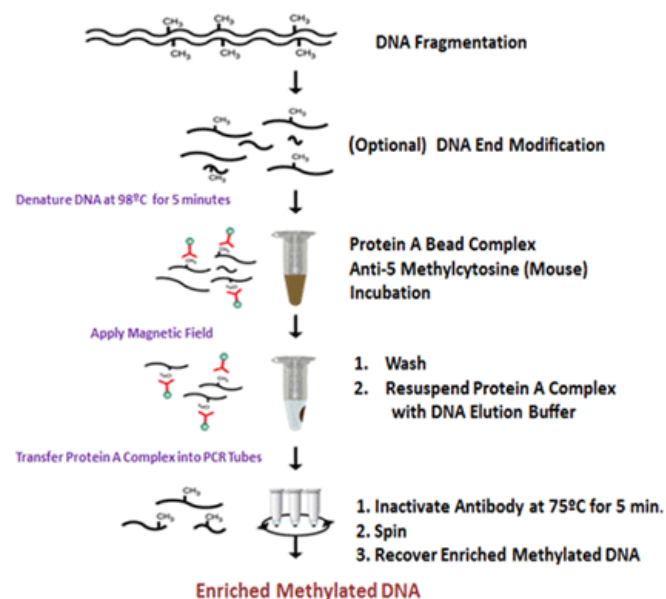


Figure demonstrates efficient enrichment of methylated vs. non methylated DNA by using the BIOARRAY™DNA IP kit. Salmon sperm genomic DNA “spiked” with non methylated/methylated control DNA was processed with the kit. The eluted DNA was amplified by PCR using the supplied control primers. Digestion of the amplicons with Nco I produced two 175bp fragments for the methylated DNA control or one 350bp fragment for the non methylated DNA control. The results show an efficient enrichment of methylated versus non methylated DNA by the kit.

Handling & Storage

| | |
|-------------------|---|
| Use/Stability | With proper storage, good for 12 months upon receipt. |
| Handling | Store each reagent based on recommended temperatures. |
| Long Term Storage | -20°C |
| Shipping | Dry Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|----------|---|
| Contents | Protein A Beads, Anti-5 methylcytosine, Control DNA, Control primer 1 , Control primer 2, IP Buffer, DNA denaturing Buffer, Control mouse DNA (methylated/nonmethylated), DNA elution buffer, Rods (magnetic) |
|----------|---|



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)