

Express DNA methylation kit

**Fastest method of complete bisulfite
conversion of DNA for methylation treatment**

The Express DNA methylation kit can be used for rapid and reliable bisulfite treatment and conversion of DNA for methylation analysis. Simply add the Express Conversion Reagent for the reaction to occur. With the addition of heat, DNA denaturation is facilitated. Desulphonation and clean-up of the converted DNA is performed using a unique low-elution spin column. High yield, converted DNA is ideal for PCR, array, bisulfite and next generation sequencing, etc.

Cytosine methylation is a naturally occurring base modification, in both prokaryotic and eukaryotic organisms. It involves the addition of a methyl group to the fifth carbon position of the cytosine pyrimidine ring via a methyltransferase enzyme. It has been demonstrated that aberrant DNA methylation is a widespread phenomenon in cancer and may be among the earliest changes to occur during oncogenesis. DNA methylation has also been shown to play a central role in gene imprinting, embryonic development, X-chromosome gene silencing, and cell cycle regulation.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-45001-0050	50 tests
ENZ-45001-0200	200 tests

Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use conversion reagent is added directly to DNA
- Easy to follow instructions
- High-yield
- DNA is ideal for PCR, MSP, array, bisulfite and Next-Gen sequencing

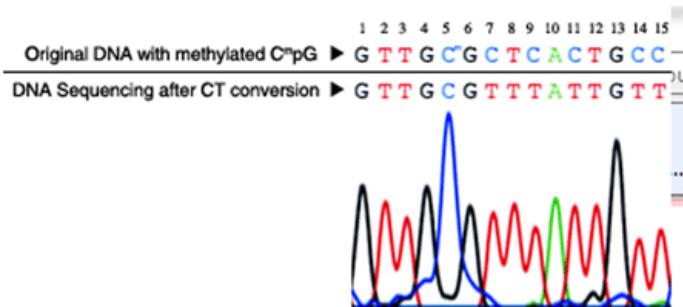


Figure: Overview of the bisulfite conversion principle. DNA template following bisulfite treatment is sequenced. Results show that methylated cytosines (position 5) remain intact while unmethylated cytosines (positions 7,11,14,15) are converted to uracil following bisulfite treatment.

Handling & Storage

Use/Stability With proper storage, good for one year upon receipt.

Handling Avoid exposure to light.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Contents Express conversion reagent, Methylation binding buffer, Methylation wash buffer concentrate, Methylation elution buffer , IC column, IC collection tube, L-Desulphonation reagent

Last modified: May 29, 2024



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)