

Express DNA methylation kit (deep well)

Fastest method for complete, high-throughput (96-well) bisulfite conversion of DNA for methylation analysis.

The Express DNA methylation kit (deep well) features high-throughput (96-well) bisulfite treatment and conversion of DNA for methylation analysis. No preparation is necessary when using the Express Conversion Reagent. Simply add this reagent to a DNA sample and let the reaction proceed to completion. DNA denaturation and bisulfite conversion processes are combined with added heat to facilitate rapid denaturation. Desulphonation and clean-up of the converted DNA is performed using a unique 96-well spin-plate. High yield, converted DNA is ideal for PCR, array, bisulfite and next generation sequencing.

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.

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

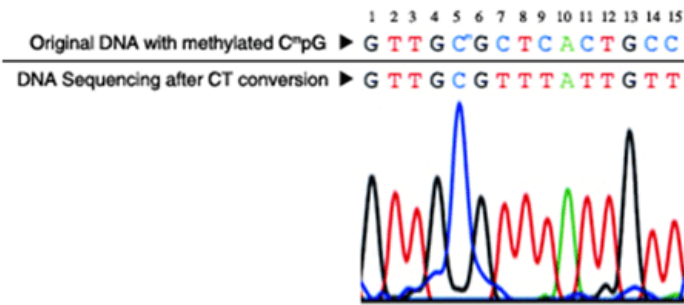


Figure: Overview of the bisulfite conversion principle. DNA template following bisulfite treatment is sequenced. Results show that methylated cytosines (position 5) remain intact while unmethylated cytosines (positions 7,11,14,15) are converted to uracil following bisulfite treatment.

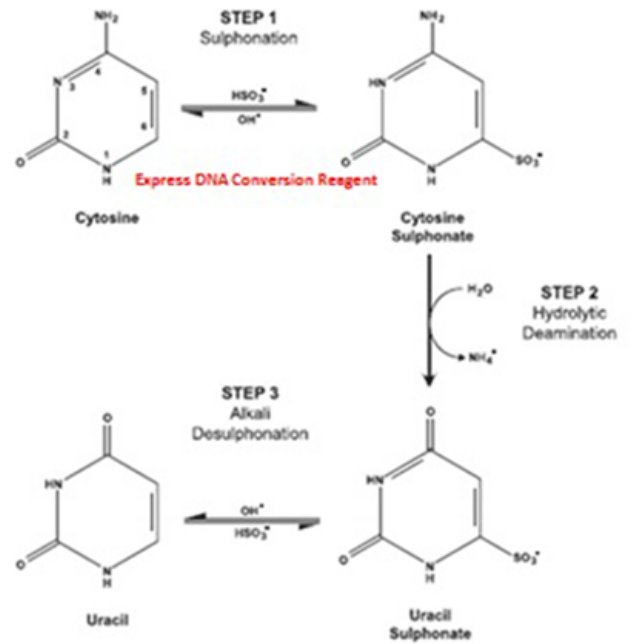


Figure: Overview of the bisulfite conversion process. Steps 1 and 2 (bisulfite conversion) occur in just one step when using the Express methylation kit. Step 3 occurs when the DNA is bound to the column matrix. DNA needs to be denatured for the reaction to be completely processed.

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, L-Desulphonation reagent, Methylation elution buffer, Express binding plate (deep), Express conversion plates, Collection plates, Elution plates
-----------------	---

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)