

MethylDetector™

simplified bisulfite conversion of DNA with easily verified results

Active Motif's MethylDetector™ Bisulfite Modification Kit simplifies analysis of DNA methylation. It comes complete with optimized reagents for performing DNA conversion with bisulfite, plus time-saving DNA purification columns and positive control PCR primers to validate your results.

DNA methylation is a naturally occurring event that affects cell function by altering gene expression. Many DNA methylation analysis methods begin by using bisulfite to convert unmethylated cytosines to uracils. However, bisulfite conversion can be technically challenging, and it is desirable to confirm that the process was successful before spending time and money on sample analysis. To help ensure your success, the MethylDetector Kit provides optimized conversion reagents, an easy-to-use protocol and positive control PCR primers that are specific for bisulfite-converted DNA.

Why use MethylDetector?

- Works efficiently with high G/C content sequences and uncut DNA
- Reproducible assay consistently provides 99% conversion efficiency of unmethylated cytosines
- Optimized reagents and protocol with proven positive controls
- Combined thermal denaturation and conversion reaction eliminates NaOH-mediated denaturation and streamlines procedure
- DNA purification columns eliminate the need for separate precipitation and desulfonation steps
- High yield of converted DNA is ideal for downstream analysis

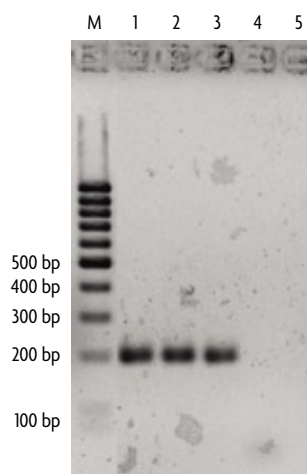


Figure 1: Agarose gel analysis of PCR products generated with MethylDetector.

Three different DNA conversions were performed (Lanes: 1-3) and compared to an unconverted DNA control (Lane: 5) and to a no DNA control (Lane: 4). The presence of PCR product in only the converted samples demonstrates the conversion efficiency and reproducibility of the MethylDetector Kit.

Product	Format	Catalog No.
MethylDetector™	50 rxns	55001

CONTENTS

Conversion Reagent, Denaturation Reagent, Hydroquinone, DNA purification columns and collection tubes, DNA Binding, Wash and Elution Buffers, Positive control PCR primers and 10X PCR Buffer.