

Methylated DNA Immunoprecipitation (MeDIP) Kit - DNA ab117133

11 References 画像数 3

製品の概要

製品名	Methylated DNA Immunoprecipitation (MeDIP) Kit - DNA
全工程の試験時間	3h 00m
製品の概要	<p>Methylated DNA Immunoprecipitation (MeDIP) Kit - DNA (ab117133) enables the user to enrich methylated DNA by using an antibody specific to methyl cytosine (5-mC) to immunoprecipitate methylated genomic DNA (<i>MeDIP</i>). The enriched methylated fractions can then be used in various downstream applications including qualitative and quantitative PCR, as well as southern blot and DNA microarray.</p> <p>DNA methylation is a covalent modification of the cytosine ring at the 5' position of a CpG dinucleotide which leads to epigenetic inactivation of genes when found in 5'-CpG-3'dinucleotides within promoters or in the first exon of genes. It has been demonstrated that DNA methylation plays an important role in the regulation of gene expression, tumorigenesis, and other genetic and epigenetic diseases, and that alterations in DNA methylation patterns are associated with cancer.</p>
特記事項	<p>NOTE: the size of this kit is based on number of TESTS not SAMPLES. A test simply refers to a single assay well. If you are not sure which size is most suitable for you, contact our Scientific Support Team.</p>

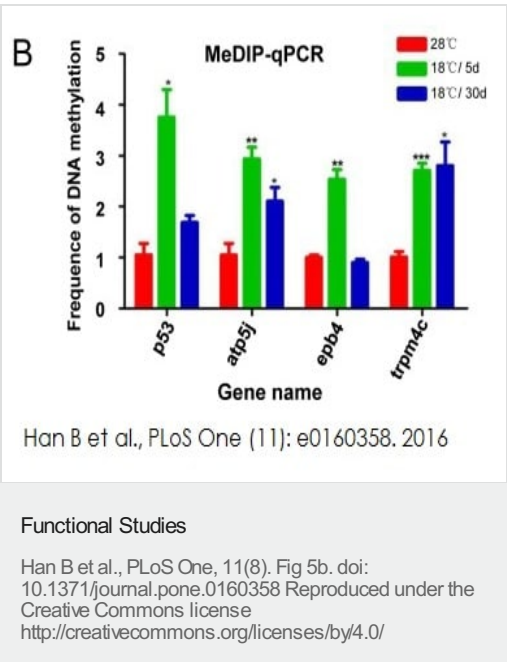
製品の特性

保存方法 Store at +4°C. Please refer to protocols.

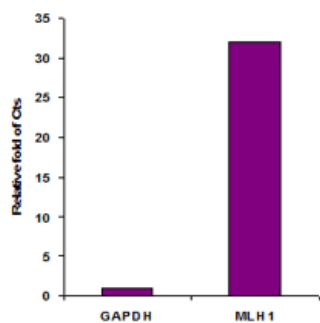
内容	24 tests	48 tests
8-Well Assay Strips (with Frame)	3 units	6 units
8-Well Strip Caps	3 units	6 units
Anti-5-Methylcytosine (1 mg/mL)	1 x 25µl	1 x 50µl
Antibody Buffer	1 x 8ml	1 x 16ml
Binding Buffer	1 x 5ml	1 x 8ml

内容	24 tests	48 tests
DNA Release Buffer	1 x 2ml	1 x 4ml
Elution Buffer	1 x 0.6ml	1 x 1.2ml
F-Collection Tube	30 units	50 units
F-Spin Column	30 units	50 units
Normal Mouse IgG (1 mg/mL)	1 x 10μl	1 x 20μl
Proteinase K (10 mg/mL)	1 x 25μl	1 x 50μl
Reaction Buffer	1 x 4ml	1 x 8ml
Wash Buffer	1 x 16ml	2 x 16ml

画像



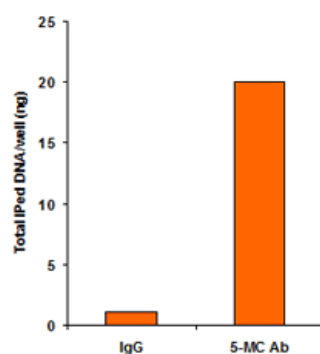
MeDIP was performed using 1μg of DNA with ab117133. Real-time PCR was carried out with MeDIP DNAs and the frequency of DNA methylation of immunoprecipitated DNA vs input DNA was calculated.



Captured methylated DNA was used for analyzing methylation level of GAPDH and MLH1 promoter with the use of primers and probes specific to GAPDH and MLH1 promoters, respectively.

Functional Studies - Methylated DNA

Immunoprecipitation (MeDIP) Kit - DNA (ab117133)



For enrichment of methylated DNA using ab117133, DNA (0.5 ug) isolated from MCF-7 cells was added into the microwell. Methylated DNA was captured by 5-mC antibody prebound to the microwells.

Functional Studies - Methylated DNA

Immunoprecipitation (MeDIP) Kit - DNA (ab117133)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors