


Methylated DNA Immunoprecipitation (MeDIP) ChIP Kit

ab117135

2 References

製品の概要

製品名	Methylated DNA Immunoprecipitation (MeDIP) ChIP Kit
サンプルの種類	Adherent cells, Suspension cells
全工程の試験時間	4h 00m
種交差性	交差種: Mouse, Human 交差が予測される動物種: Rat 

製品の概要	<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> <p>Methylated DNA Immunoprecipitation (MeDIP) ChIP Kit (ab117135) enables the user to enrich methylated DNA directly from cells 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>
-------	--

製品の特性

保存方法	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 15ml	1 x 30ml
Binding Buffer	1 x 5ml	1 x 8ml

内容	24 tests	48 tests
ChIP Dilution Buffer	1 x 4ml	1 x 6ml
DNA Release Buffer	1 x 2ml	2 x 2ml
Elution Buffer	1 x 0.6ml	1 x 1.2ml
F-Collection Tube	30 units	50 units
F-Spin Column	30 units	50 units
Lysis Buffer	1 x 4ml	1 x 6ml
Normal Mouse IgG (1 mg/mL)	1 x 10µl	1 x 15µl
Pre-Lysis Buffer	1 x 4ml	1 x 8ml
Proteinase K (10 mg/mL)	1 x 25µl	1 x 50µl
Reverse Buffer	1 x 2ml	2 x 2ml
Wash Buffer	1 x 28ml	2 x 28ml

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