


Bisulfite Conversion Kit - Whole Cell ab117126

画像数 1

製品の概要

製品名	Bisulfite Conversion Kit - Whole Cell
サンプルの種類	Saliva, Urine, Tissue, Adherent cells, Suspension cells
全工程の試験時間	3h 00m
種交差性	交差種: Human 交差が予測される動物種: Mammals 

製品の概要

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 is well demonstrated that DNA methylation plays an important role in the regulation of gene expression, tumorigenesis, and other genetic and epigenetic diseases.

Most methods for DNA methylation detection require a prior bisulfite-based DNA modification. *Bisulfite-based DNA modification* is used to discriminate between cytosine and methylated cytosine, in which DNA is treated with bisulfite salt to convert cytosine residues to uracil in a single-stranded DNA, while methylated cytosine remains the same.

Bisulfite Modification Kit - Whole Cell (ab117126) allows the user to perform DNA methylation analysis and modify DNA directly from **cells** or **tissues** within only 3 hours. ab117126 is specifically designed for DNA methylation research using minute amounts of starting materials including cells cultured in 96-well/384-well plates, tissue section samples, microdissection samples, tissue biopsy and early embryonic cells/oocytes.

Modified DNA is eluted and suitable for real time MS-PCR as well as for all techniques currently used for the analysis of DNA methylation; including conventional MS-PCR, bisulfite sequencing, pyrosequencing, and methylation microarray.

Not sure if this is the right kit for you? Check out our [bisulfite modification guide](#) for more information.

特記事項

法規制情報

医薬用外劇物

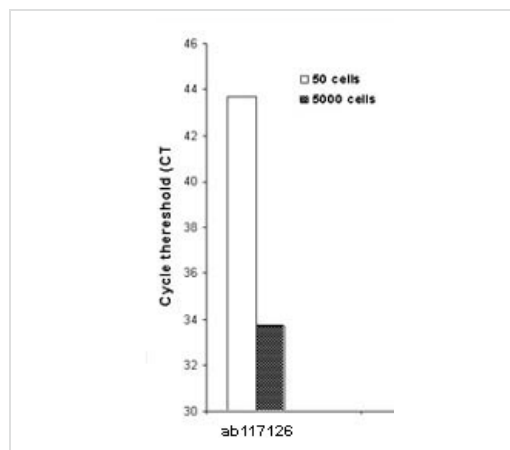
製品の特性

保存方法

Please refer to protocols.

内容	40 tests	80 tests
Balance Buffer	1 x 0.4ml	1 x 0.8ml
Cell Collection Buffer	1 x 1ml	1 x 2ml
Digestion Powder	1 vial	1 vial
Digestion Solution	1 x 0.1ml	1 x 0.2ml
DNA Binding Buffer	1 x 14ml	1 x 28ml
DNA Modification Buffer	1 x 5ml	1 x 10ml
DNA Modification Powder	4 vials	8 vials
F-Collection Tube	1 x 40 units	1 x 80 units
F-Spin Column	1 x 40 units	1 x 80 units
Modified DNA Elution	1 x 1ml	1 x 2ml

画像



Functional Studies - Bisulfite Modification Kit -
Whole Cell (ab117126)

MCF-7 cells were modified using ab117126. 10 μ l of modified DNA were eluted and 2 μ l of elution were used in real time PCR. A pair of primers and a probe designed to amplify both methylated and unmethylated alleles of β -actin was used.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Promise 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