

# Methylated Control DNA Kit ab117125

## 製品の概要

**製品名** Methylated Control DNA Kit

**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.

A methylation-positive control is a must tool for a successful performance of gene methylation studies.

Abcam's Methylated Control DNA Kit (ab117125) provides a methylation-positive control for methylation studies. The kit includes the enzymatically methylated human genomic DNA and all components for DNA modification. Methylated human DNA needs to be modified before it is used as a positive control in MS-PCR. DNA modification reagents included in the kit are sufficient for modifying 5 µg of methylated human genomic DNA (10 modifications).

This product is intended for use with the Abcam's DNA modification kit series. It can also be used with other DNA modification or methylation kits.

**アプリケーション** **適用あり:** Functional Studies

## 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	10 tests	20 tests
Balance Solution	1 x 0.08ml	1 x 0.15ml
DNA Binding Buffer	1 x 4ml	1 x 8ml
DNA Modification Solution	1 x 1.5ml	1 x 3ml
DNA Modifier Powder	1 x 0.7g	1 x 1.4g

内容	10 tests	20 tests
Elution Buffer	1 x 0.5ml	1 x 1ml
F-Collection Tube	10 units	20 units
F-Spin Column	10 units	20 units
Methylated Human Genomic DNA (0.1 mg/mL)	1 x 5µg	1 x 10µg

## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab117125** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

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

## 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 <http://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