

# Post-Bisulfite DNA Library Preparation Kit (For Illumina®) ab185906

## 1 References [画像数 1](#)

### 製品の概要

---

製品名	Post-Bisulfite DNA Library Preparation Kit (For Illumina®)
検出感度	1 ng
検出範囲	0.5 ng - 1000 ng
全工程の試験時間	5h 00m
製品の概要	ab185906 is a complete set of optimized reagents to prepare a DNA library - after successful bisulfite conversion - for various Illumina® platform-based bisulfite sequencing (bisulfite-seq) assays, such as whole genome bisulfite sequencing (WGBS), oxidative bisulfite sequencing (oxBS-seq), reduced representative bisulfite sequencing (RRBS), and other bisulfite-based next generation sequencing applications. The optimized protocol and components of the kit allow both non-barcoded (singleplexed) and barcoded (multiplexed) DNA libraries to be quickly constructed using sub-nanogram input concentrations of DNA since the DNA is first bisulfite-converted and then used for library preparation.

Illumina® is a registered trademark of Illumina, Inc.

### 特記事項

Several methods, including whole genome bisulfite sequencing (WGBS) and reduced representation bisulfite sequencing (RRBS), are currently used for genome-wide DNA methylation analysis. These methods convert unmethylated cytosines to uracil while 5-methylcytosines remain unchanged by the bisulfite treatment. This allows epigenetic differences to become genetic differences, which can be subsequently detected via sequencing at the single-based resolution level and on a genome-wide scale. However, practical use with such current methods are not ideal as they (1) need large amounts of DNA (>1 µg) as input material, which is difficult to prepare from limited biological samples such as tumor biopsy, early embryos, embryonic tissues, and circulating DNA; (2) require DNA to first be sheared and then ligated to adapters, followed by bisulfite conversion (post-ligation bisulfite conversion), which causes substantial amounts of DNA fragments contained in the adapter-DNA fragment constructs to be broken and thereby forms mono-tagged templates that become removed during library enrichment; and (3) are time-consuming (2 days). The Post-Bisulfite DNA Library Preparation Kit (For Illumina®) is designed to overcome these weaknesses.

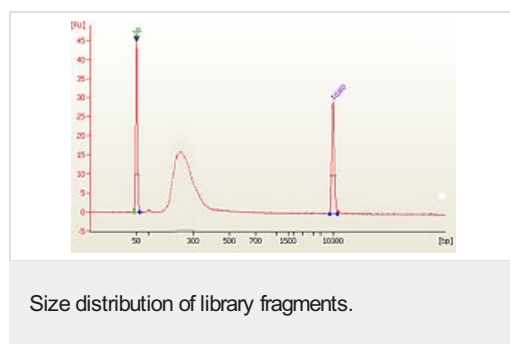
## 製品の特性

### 保存方法

Please refer to protocols.

内容	12 tests	24 tests
10X dA-Tailing Buffer	1 x 40 $\mu$ l	1 x 80 $\mu$ l
10X End Repair Buffer	1 x 40 $\mu$ l	1 x 80 $\mu$ l
2X HiFi PCR Master Mix	1 x 160 $\mu$ l	1 x 320 $\mu$ l
2X Ligation Buffer	1 x 250 $\mu$ l	1 x 500 $\mu$ l
5X Conversion Buffer	1 x 50 $\mu$ l	1 x 100 $\mu$ l
Adaptors (50 $\mu$ M)	1 x 15 $\mu$ l	1 x 30 $\mu$ l
Conversion Enzyme Mix	1 x 15 $\mu$ l	1 x 30 $\mu$ l
Conversion Primer	1 x 26 $\mu$ l	1 x 52 $\mu$ l
Elution Buffer	1 x 1ml	1 x 2ml
End Repair Enzyme Mix	1 x 25 $\mu$ l	1 x 50 $\mu$ l
Klenow Fragment (3'-5' exo-)	1 x 15 $\mu$ l	1 x 30 $\mu$ l
MQ Binding Beads	1 x 1.8ml	1 x 3.6ml
Primer I (10 $\mu$ M)	1 x 15 $\mu$ l	1 x 30 $\mu$ l
Primer U (10 $\mu$ M)	1 x 15 $\mu$ l	1 x 30 $\mu$ l
T4 DNA Ligase	1 x 15 $\mu$ l	1 x 30 $\mu$ l

### 画像



Size distribution of library fragments as demonstrated by a post-bisulfite DNA library constructed using ab185906 Post-Bisulfite DNA Library Preparation Kit (For Illumina®) from 10 ng of input DNA.

**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