

### Bisulfite-Seq High Sensitivity Kit (For Illumina®) ab185907

画像数 1

医薬用外劇物

#### 製品の概要

|          |   |
|----------|---|
| 製品名      | Bisulfite-Seq High Sensitivity Kit (For Illumina®)  |
| 検出感度     | < 0.5 ng  |
| 全工程の試験時間 | 8h 00m  |
| 製品の概要    | ab185907 is designed to carry out bisulfite conversion, followed by a "post-bisulfite" library preparation process for Illumina® platform-based bisulfite sequencing, all in one kit. The DNA library is constructed directly from bisulfite-converted DNA generated from a small amount of input DNA (500 pg to 500 ng). Intended applications include whole genome bisulfite sequencing, oxidative bisulfite sequencing, reduced representative bisulfite sequencing, and various other bisulfite-next generation sequencing techniques. The optimized protocol and components of the kit allow the DNA to be bisulfite converted and fragmented simultaneously followed by quick non-barcoded (singleplexed) and barcoded (multiplexed) library construction using sub-nanogram quantities of bisulfite converted DNA. |

#### 特記事項

DNA methylation occurs by the covalent addition of a methyl group (CH<sub>3</sub>) at the 5-carbon of the cytosine ring, resulting in 5-methylcytosine (5-mC). DNA methylation is essential in regulating gene expression in nearly all biological processes including development, growth, and differentiation. Alterations in DNA methylation have been demonstrated to cause a change in gene expression. Genome-wide analysis of DNA methylation could provide valuable information for discovering epigenetic markers used for disease diagnosis, and potential targets used for therapeutics. Bisulfite sequencing via next-generation sequencing technologies allow for high volume, lower cost output of DNA sequence data towards a better understanding of DNA methylation.

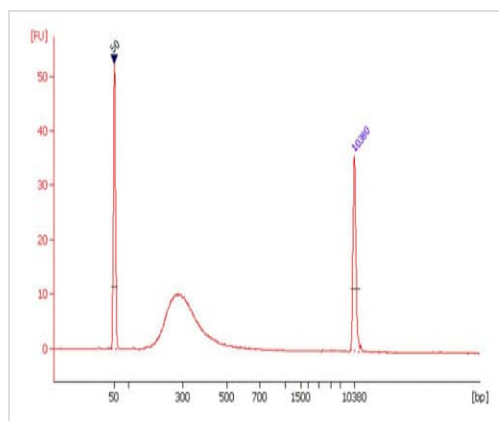
#### 製品の特性

保存方法 Please refer to protocols.

| 内容                    | 24 tests | 12 tests |
|-----------------------|----------|----------|
| 10X dA-Tailing Buffer | 1 x 80µl | 1 x 40µl |
| 10X End Repair Buffer | 1 x 80µl | 1 x 40µl |

| 内容                           | 24 tests     | 12 tests     |
|------------------------------|--------------|--------------|
| 2X HiFi PCR Master Mix       | 1 x 320µl    | 1 x 160µl    |
| 2X Ligation Buffer           | 1 x 500µl    | 1 x 250µl    |
| 5X Conversion Buffer         | 1 x 100µl    | 1 x 50µl     |
| Adaptors (50 µM)             | 1 x 30µl     | 1 x 15µl     |
| Conversion Enzyme Mix        | 1 x 30µl     | 1 x 15µl     |
| Conversion Primer            | 1 x 52µl     | 1 x 26µl     |
| Desulphonation Solution      | 1 x 140µl    | 1 x 70µl     |
| DNA Binding Solution         | 1 x 12ml     | 1 x 6ml      |
| Elution Buffer               | 1 x 2ml      | 1 x 1ml      |
| Elution Solution             | 1 x 1ml      | 1 x 0.5ml    |
| End Repair Enzyme Mix        | 1 x 50µl     | 1 x 25µl     |
| F-Collection Tube            | 1 x 30 units | 1 x 15 units |
| F-Spin Column                | 1 x 30 units | 1 x 15 units |
| Klenow Fragment (3'-5' exo-) | 1 x 30µl     | 1 x 15µl     |
| Modification Buffer          | 1 x 6ml      | 1 x 3ml      |
| Modification Powder          | 1 x 4 vials  | 1 x 2 vials  |
| MQ Binding Beads             | 1 x 3.6ml    | 1 x 1.8ml    |
| Primer I (10 µM)             | 1 x 30µl     | 1 x 15µl     |
| Primer U (10 µM)             | 1 x 30µl     | 1 x 15µl     |
| T4 DNA Ligase                | 1 x 30µl     | 1 x 15µl     |

画像



Size distribution of library fragments. Post-bisulfite DNA library was prepared from 10 ng of input DNA using ab185907.

Size distribution of library fragments

**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