

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

1 图像

概述

| | |
|------|---|
| 产品名称 | Bisulfite-Seq High Sensitivity试剂盒(For Illumina®) |
| 灵敏度 | < 0.5 ng |
| 检测时间 | 8h 00m |
| 产品概述 | 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. |

说明

DNA methylation occurs by the covalent addition of a methyl group (CH₃) 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.

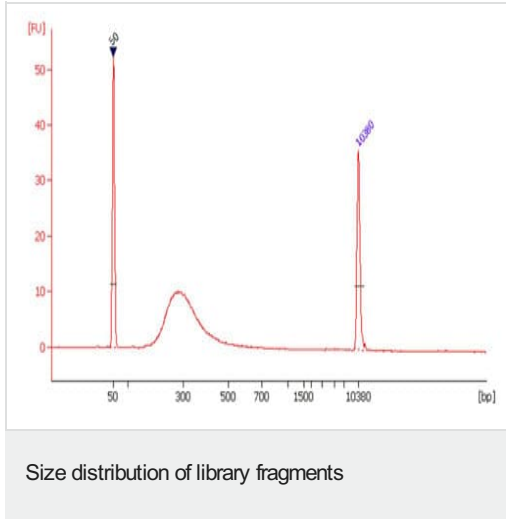
性能

存放说明 Please refer to protocols.

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

| 组件 | 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 |

图片



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

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors