

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

[1 References](#) [1 Image](#)

Overview

Product name	Post-Bisulfite DNA Library Preparation Kit (For Illumina®)
Sensitivity	1 ng
Range	0.5 ng - 1000 ng
Assay time	5h 00m
Product overview	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 representation 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.

Notes	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.
--------------	--

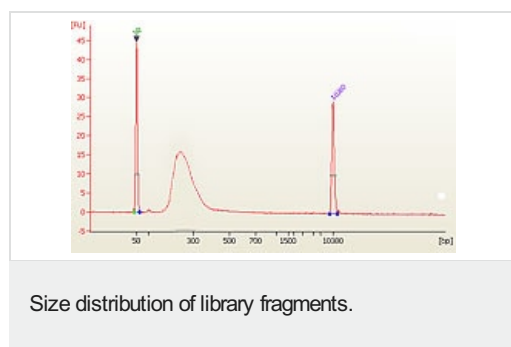
Properties

Storage instructions

Please refer to protocols.

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

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

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