AMPINEXT™ High-Sensitivity Bisulfite-Seq Kit (Illumina)

Designed to bisulfite-convert DNA and prepare an Illumina-based library for bisulfite

The AMPINETY To Maigh Sensitivity Bisulfite-Seq Kit (Illumina) is designed to bisulfite-convert DNA and prepare an Illumina-based library for bisulfite sequencing, all in one kit. Intended applications include whole genome (WGBS), oxidative (oxBS-seq), reduced representation (RRBS), and other bisulfite-next generation sequencing. The optimized protocol and components of the kit allow subnanogram amounts of DNA to be bisulfite converted and fragmented simultaneously followed by quick non-barcoded (singleplexed) and barcoded (multiplexed) library construction in less than 7 hours.

Ordering Information

Order Online »

ENZ-GEN508-0012	12Reactions
ENZ-GEN508-0024	24Reactions

Manuals, SDS & CofA

View Online »

- Innovative method Allows for simultaneous bisulfite conversion and size-appropriate DNA fragmentation.
- Fast and streamlined procedure
 - Ready-to-use library DNA within 6 hours and 30 minutes.
- Complete conversion Converts unmethylated cytosine into uracil (>99%).
- High sensitivity and efficiency -Innovative adaptor ligation of bisulfite DNA eliminates loss of fragments and selection bias.
- Extremely convenient The kit contains all required components for each step of DNA library preparation,
- Minimal bias Ultra HiFi amplification.
- Broad sample suitability Starting materials can be genomic
 DNA isolated from various
 tissue/cell samples.

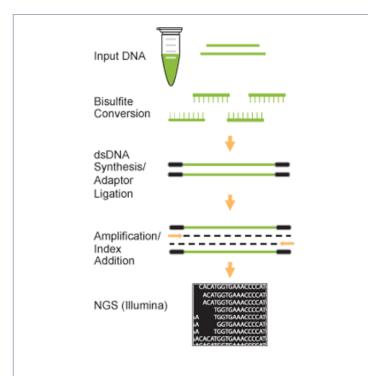


Figure 1. Schematic Procedure for using the AMPINEXT High-Sensitivity Bisulfite-Seq Kit (Illumina)

Handling & Storage

Use/Stability Upon receipt: (1) Store the following components at -20°C immediately: 5X Reaction

Buffer, Reaction Enzyme Mix, Adaptor-F, Adaptor-S, 2X HiFi PCR Master Mix, Primer U, Primer I, and Elution Buffer. (2) Store the following components at 4°C: MQ Binding

Beads. (3) Store all other components at room temperature away from light.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application NotesDesigned to bisulfite-convert DNA and prepare an

Illumina-based library for bisulfite sequencing, all in one

kit.

Assay Time <7 hours

Contents Modification Buffer

Modification Powder
DNA Binding Solution

Desulphonation Solution

Elution Solution F-Spin Column

F-Collection Tube

5X Reaction Buffer Reaction Enzyme Mix

Adaptor-F (10 μM) Adaptor-S (10 μM)

MQ Binding Beads

2X HiFi PCR Master Mix

Primer U (10 µM) Primer I (10 µM) Elution Buffer

Last modified: May 29, 2024

