

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

Designed to bisulfite-convert DNA and prepare an Illumina-based library for bisulfite sequencing, all in one kit

The AMPINEXT™ High-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.

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.

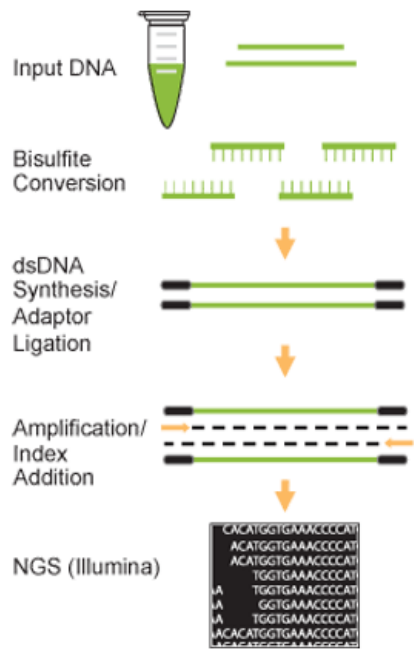


Figure 1. Schematic Procedure for using the AMPINEXT High-Sensitivity Bisulfite-Seq Kit (Illumina)

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 Notes Designed 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com