AMPINEXT™ Post-Bisulfite DNA Library Preparation Kit (Illumina)

The AMPINEXT'S Post Bis little DNA library Greparation Kit is a complete set of optimized reagents to prepare a DNA library - after successful bisulfite conversion -- for various Illumina platform-based bisulfite sequencing (bisulfite set) assays, such as whole genome bisulfite pagueroing (WSBS) exidative bisulfite sequencing (oxBs-seq), reduced representation bisulfite sequencing (RRBS), and other bisulfite-based next (DISUTITE-SEQ) assays, 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 subnanogram input concentrations of DNA since the DNA is first bisulfite-converted and then used for library preparation.

Ordering Information

Order Online »

ENZ-GEN512-0012	12Reactions
ENZ-GEN512-0024	24Reactions

Manuals, SDS & CofA

View Online »

- Allows bisulfite-converted DNA to be used directly for ligation.
- Fast 5-hour procedure, from input starting material to library amplification.
- Gel-free size selection/purification
- High sensitivty and efficiency -direct ligation of adapter to bisulfite-converted DNA fragments.
- Comprehensive set of components to accommodate each step of DNA library preparation.
- Ultra HiFi amplification enables achievement of reproducibly high yields of DNA libraries.

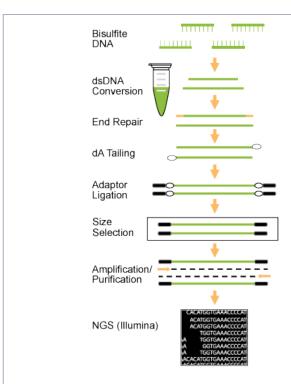


Figure 1. Schematic Procedure for using the AMPINEXT Post-Bisulfite DNA Library Preparation Kit (Illumina) Schematic

Handling & Storage

Use/Stability Upon receipt, store the following components at -20°C immediately: 5X Conversion

Buffer, Conversion Enzyme Mix, Conversion Primer, 10X End Repair Buffer, End Repair Enzyme Mix,10X dA-Tailing Buffer, Klenow Fragment (3'-5' exo-), 2X Ligation Buffer, T4 DNA Ligase, Adaptors, 2X HiFi PCR Master Mix, Primer U, Primer I, and Elution Buffer. Store the following components at 4°C: MQ Binding Beads. Store all other components

at room temperature.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes 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.

Assay Time 5 hours

Contents 5X Conversion Buffer

Conversion Enzyme Mix

Conversion Primer

10X End Repair Buffer

End Repair Enzyme Mix

10X dA-Tailing Buffer

Klenow Fragment (3'-5' exo⁻)

2X Ligation Buffer

T4 DNA Ligase

Adaptors (50 µM)

MQ Binding Beads

2X HiFi PCR Master Mix

Primer U (10 µM)

Primer I (10 µM)

Elution Buffer

Last modified: February 7, 2025



usa@enzolifesciences.com