

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

The AMPINEXT™ Post-Bisulfite DNA Library Preparation 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-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.

Ordering Information

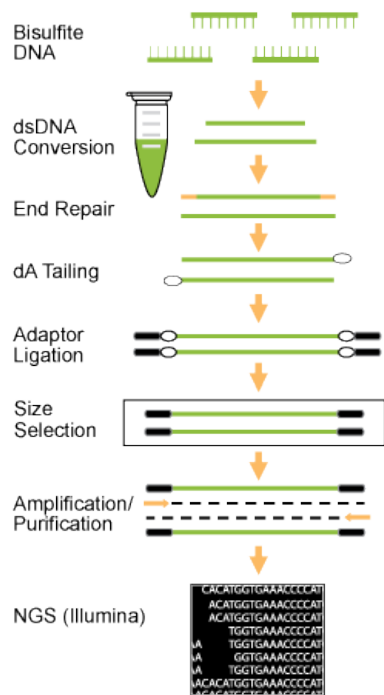
[Order Online »](#)

ENZ-GEN512-0012	12Reactions
-----------------	-------------

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 sensitivity 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.



```

CACATGGTGAACCCCAT
ACATGGTGAACCCCAT
ACATGGTGAACCCCAT
TGGTGAACCCCAT
AA TGGTGAACCCCAT
AA GGTGAACCCCAT
AA TGGTGAACCCCAT
AACACATGGTGAACCCCAT
AAACACATGGTGAACCCCAT

```

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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