

AMPINEXT™ DNA Library Preparation Kit (Illumina)

Complete set of optimized reagents to carry out a successful DNA library preparation.

The AMPINEXT™ DNA Library Preparation Kit (Illumina) is a complete set of optimized reagents to carry out a successful DNA library preparation. The kit is suitable for preparing a DNA library for next generation sequencing applications using an Illumina sequencer, which includes genomic DNA-seq, ChIP-seq, MeDIP/hMeDIP-seq, bisulfite-seq, and targeted re-sequencing. The optimized protocol and components of the kit allow both non-barcoded (singleplexed) and barcoded (multiplexed) DNA libraries to be constructed quickly with reduced bias.

Ordering Information

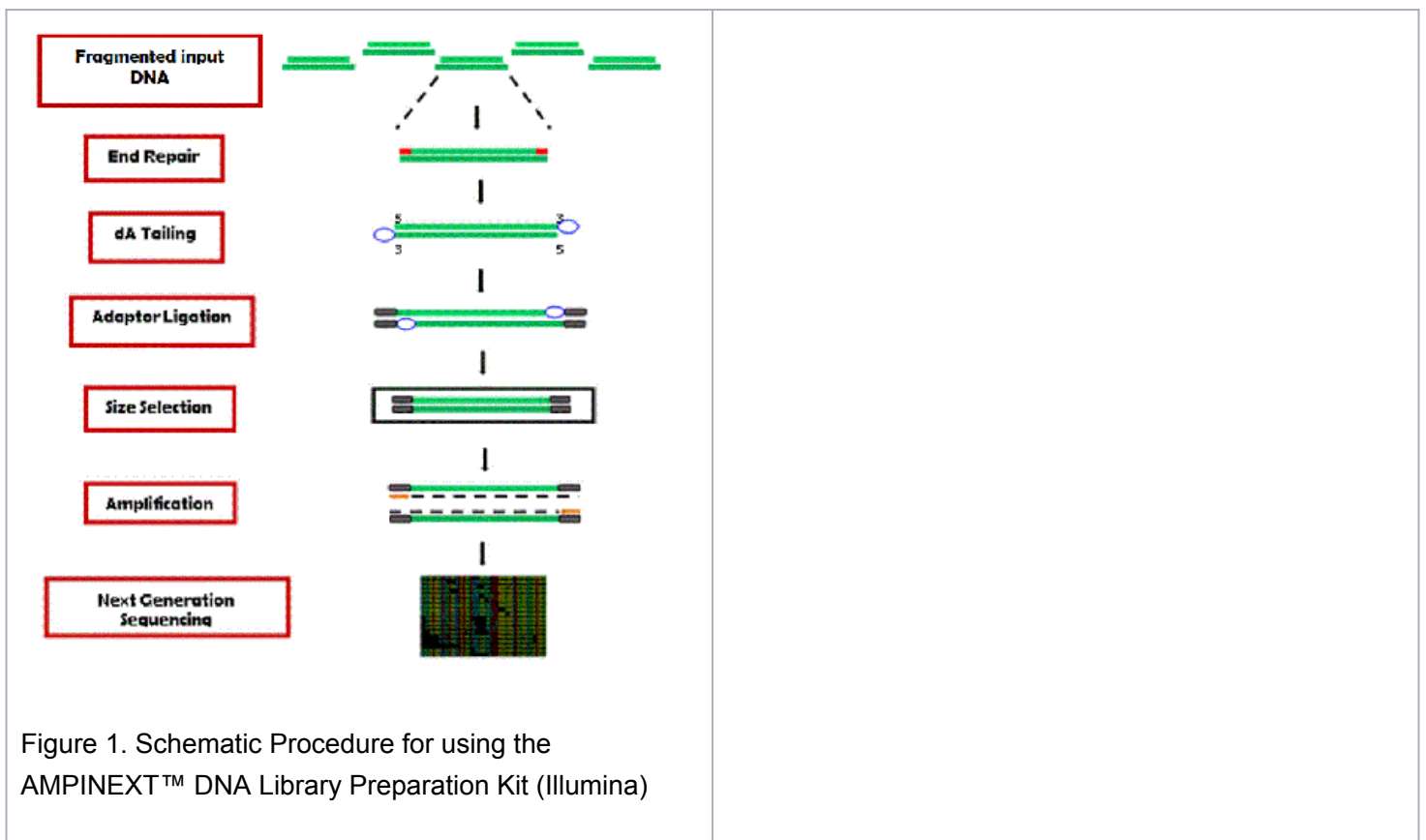
[Order Online »](#)

ENZ-GEN504-0012	12Reactions
ENZ-GEN504-0024	24Reactions

Manuals, SDS & CofA

[View Online »](#)

- **Fast and streamlined procedure**
 - The procedure from fragmented DNA to size selection is less than 2 h 30 min.
- **Highly Convenient** - The kit contains all required components for each step of DNA library preparation.
- **Minimized bias** - Ultra HiFi amplification and an optional PCR-free step.
- **Flexibility** - Use for both non-barcoded (singleplexed) and barcoded (multiplexed) DNA library preparation.



Handling & Storage

Use/Stability Upon receipt: Store the following components at -20°C immediately: 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. □

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes A complete set of optimized reagents to carry out a successful DNA library preparation.

Assay Time 2 hours 30 minutes

Contents

- 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

Technical Info / Product Notes Starting materials can include fragmented dsDNA isolated from various tissue or cell samples, dsDNA enriched from ChIP reactions, MeDIP/hMeDIP reaction, or exon capture. DNA should be relatively free of RNA since large fractions of RNA will impair end repair and dA tailing, resulting in reduced ligation capabilities. Input amount of DNA can be from 5 ng to 1 µg. For optimal preparation, the input amount should be 100 ng to 200 ng. For amplification-free, 500 ng or more is needed.



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-
fr@enzolifesciences.com](mailto:info-fr@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)