



AMPINEXT™ High Sensitivity DNA Library Preparation Kit (Illumina)

Catalog #: *ENZ-GEN505*

ENZ-GEN505-0012 – for 12 reactions

ENZ-GEN505-0024 – for 24 reactions

For the latest product information, including support documentation, visit us online:

www.enzolifesciences.com

Technical Support (US): 800-942-0430
Technical Support (EU): +41 61 926 8989



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research use only. Not for use in diagnostic procedures. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON- INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

TABLE OF CONTENTS



Carefully note the handling and storage conditions of each kit component.

INTRODUCTION	2
BACKGROUND	2
MATERIALS SUPPLIED	3
ADDITIONAL MATERIALS NEEDED	4
STORAGE & STABILITY	4
SAFETY WARNINGS & PRECAUTIONS	5
PRINCIPLE OF THE ASSAY	5
PROCEDURE	6
TROUBLESHOOTING GUIDE	13
CONTACT INFORMATION	Error! Bookmark not defined.

INTRODUCTION

DNA library preparation is a critical step for next generation sequencing (NGS). For generating accurate sequencing data in NGS, the prepared library DNA should be sufficient in yield and of high quality. Also as NGS technology is continuously improving, DNA library preparation is required to be optimized accordingly. For example, most of the currently used methods are time-consuming, expensive, inconvenient, and specifically need large amounts of DNA. These reactions result in a DNA library preparation which cannot be used for biological samples with limited amounts of starting material such as tumor biopsy, early embryos, embryonic tissues and circulating DNA. In addition, the amount of DNA enriched by ChIP or MeDIP/hmeDIP is often at low or sub-nanogram levels which causes insufficient DNA library yields. To address this issue, we offer the AMPINEXT™ High Sensitivity DNA Library Preparation Kit (Illumina). This kit has the following features:

- High sensitivity and flexibility: Can be used for both non-barcoded (singleplexed) and barcoded (multiplexed) DNA library preparation. The amount of input DNA can be as low as 0.2 ng with a range from 0.2 to 100 ng. Various dsDNA can be used, which includes limited amounts of fragmented dsDNA isolated from various tissue or cell samples, dsDNA enriched from ChIP reactions, and dsDNA enriched from MeDIP/hMeDIP reactions or exon capture.
- Fast and streamlined procedure: the procedure from fragmented DNA to size selection is less than 1 h 30 min. No clean-up is required between each step and all reactions take place in the same tube, thereby saving time and preventing handling errors, as well as loss of valuable samples. Gel-free size selection further reduces the preparation time.
- The most convenient for use: the kit contains all required components for each step of DNA library preparation, which are sufficient for end polishing, ligation, clean-up, size selection and library amplification, thereby allowing the library preparation to be the most convenient with reliable and consistent results.
- Minimized bias: Ultra HiFi amplification enables to reproducibly achieve high yields of DNA library with minimal sequence bias and low error rates.

BACKGROUND

Uses: The AMPINEXT™ High Sensitive DNA Library Preparation Kit (Illumina) is suitable for preparing a DNA library using sub-nanogram amounts of DNA input for next generation sequencing applications using an Illumina sequencer. These applications include genomic DNA-seq, ChIP-seq, MeDIP/hMeDIP-seq,

classical 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.

Starting Material and Input Amount: Starting materials can include fragmented dsDNA isolated from various tissue or cell samples, dsDNA enriched from a ChIP reaction, MeDIP/hMeDIP reaction or exon capture. DNA should be relatively free of RNA because large fractions of RNA will impair end repair and dA-tailing, resulting in reduced ligation capabilities. The input amount of DNA can be from 0.2 ng to 100 ng. For optimal preparation, the input amount should be 10 ng to 50 ng.

MATERIALS SUPPLIED

Components	12 reactions (ENZ-GEN505 - 0012)	24 reactions (ENZ-GEN505- 0024)	Storage Upon Receipt
10X End Polishing Buffer*	30 µl	60 µl	-20°C
End Polishing Enzyme Mix*	13 µl	26 µl	-20°C
End Polishing Enhancer	13 µl	26 µl	-20°C
2X Ligation Buffer*	250 µl	500 µl	-20°C
T4 DNA Ligase*	15 µl	30 µl	-20°C
Adaptors (50 µM)*	15 µl	30 µl	-20°C
MQ Binding Beads*	1.6 ml	3.2 ml	4°C
2X HiFi PCR Master Mix*	160 µl	320 µl	-20°C
Primer U (10 µM)*	15 µl	30 µl	-20°C
Primer I (10 µM)*	15 µl	30 µl	-20°C
Elution Buffer*	1000 µl	2000 µl	-20°C
Product Manual	1	1	RT

* Spin the solution down to the bottom prior to use.

ADDITIONAL MATERIALS NEEDED

The following are required but not provided:

- Vortex mixer
- Sonicator or enzymes for DNA fragmentation
- Agilent® Bioanalyzer® or comparable method to assess the quality of DNA library
- Thermocycler
- Centrifuge including desktop centrifuge (up to 14,000 rpm)
- Magnetic stand (96-well format)
- Pipettes and pipette tips
- PCR tubes or plates
- 1.5 ml microcentrifuge tubes
- 80% Ethanol
- Distilled water
- DNA sample



Storage temp

STORAGE & STABILITY

The kit is shipped on frozen ice packs at 4°C.

Upon receipt: Store the following components at -20°C immediately: **10X End Polishing Buffer, End Polishing Enzyme Mix, End Polishing Enhancer, 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.



Important/ Warning

SAFETY WARNINGS & PRECAUTIONS

1. Wear appropriate personnel protective apparel. Avoid contact with clothes and exposed skin. In case of accidental skin exposure, flush with water immediately. Consult a physician if required.
2. Use a safety pipetting device for all pipetting. Never pipet by mouth.
3. Interpretation of the results is the sole responsibility of the user.

PRINCIPLE OF THE ASSAY

The AMPINEXT™ High Sensitive DNA Library Preparation Kit (Illumina) contains all reagents required at each step of workflow for carrying out successful DNA library preparation. In the library preparation, DNA is first fragmented to appropriate size (about 300 bps in peak size). The end repair/dA tailing (end polishing) of the DNA fragments are performed simultaneously. Adaptors are then ligated to both ends of the polished DNA fragments for amplification and sequencing. Ligated fragments are size selected and purified with MQ beads, which allows quick and precise size selection of DNA. Size-selected DNA fragments are amplified with high-fidelity PCR Mix that ensures maximum yields from minimum amounts of starting material and provides highly accurate amplification of library DNA with low error rates and minimum bias.

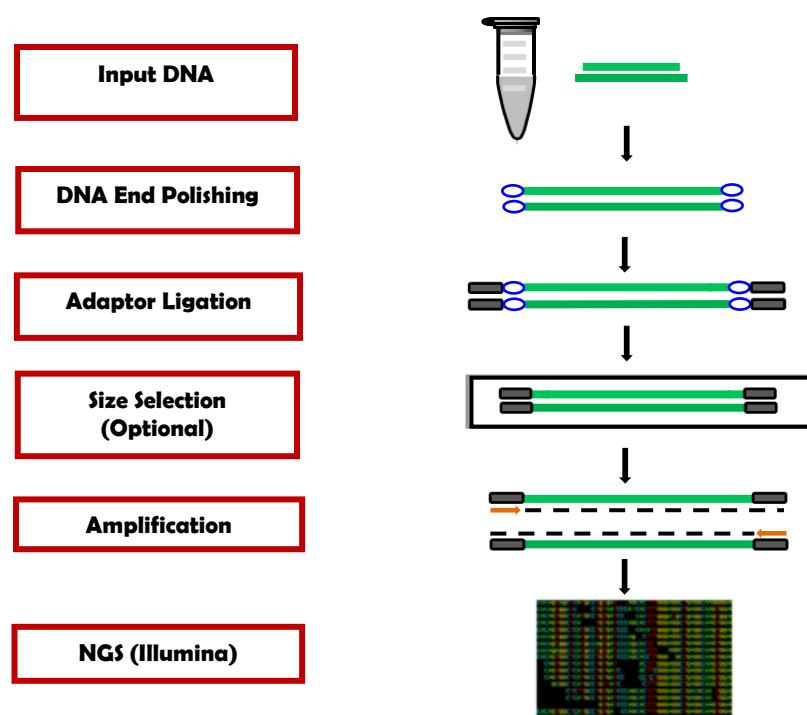


Fig 1. Workflow of the AMPINEXT™ High Sensitive DNA Library Preparation Kit

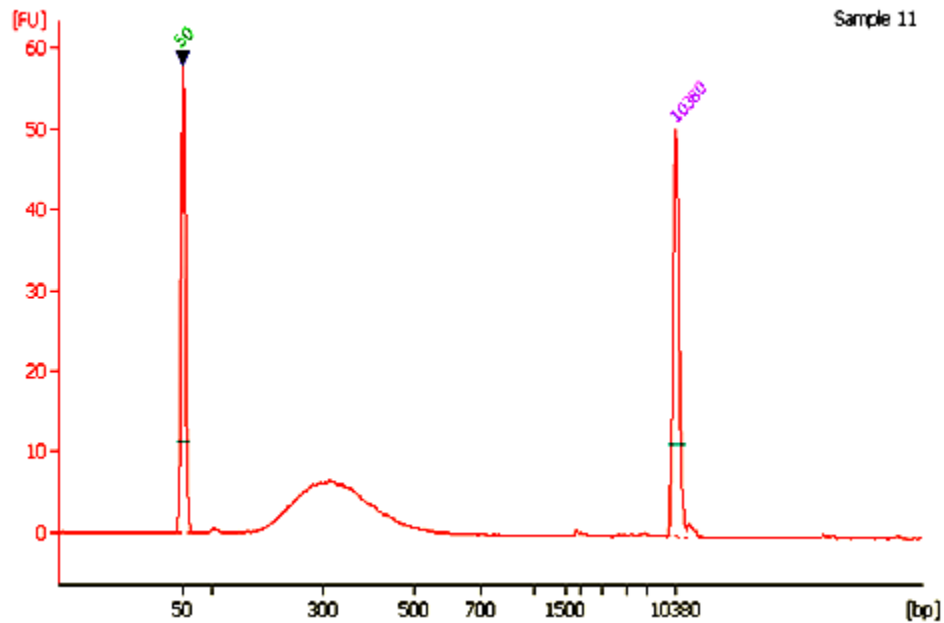


Fig 2. Size distribution of library fragments. Human placenta DNA was sheared to around 300 bps in peak size and 0.2 ng of DNA was used for DNA library preparation using AMPINEXT™ High Sensitive DNA Library Preparation Kit (Illumina)

PROCEDURE

Precautions: To avoid cross-contamination, carefully pipette the sample or solution into the tube/vials. Use aerosol-barrier pipette tips and always change pipette tips between liquid transfers. Wear gloves throughout the entire procedure. In case of contact between gloves and sample, change gloves immediately.

For the best results, please read the protocol in its entirety prior to starting your experiment.

Starting Materials

Fragmented dsDNA that is isolated from various tissues or cell samples: 0.2 ng-100 ng, optimized 10-50 ng per preparation.

dsDNA enriched from a ChIP reaction, MeDIP/hMeDIP reaction or exon capture: 0.2 ng-100 ng.

DNA should be of high quality and relatively free of RNA. RNase I can be used to remove RNA and DNA should be eluted in DNase/RNase-free water.

DNA Fragmentation

dsDNA enriched from a CHIP reaction, MeDIP/hMeDIP reaction or exon capture should already be fragmented. DNA isolated from various tissue or cell samples can be fragmented using one of the following methods. For the best results we highly recommend using a waterbath-based sonication device. The peak size of fragmented DNA should be compatible with the read length of the Illumina sequencing platform to be used. In general the peak size of fragments should be 200-300 bps.

Waterbath Sonication:

When using waterbath sonicators, please follow the supplier's instruction.

Enzymatic Shearing:

The DNA can also be sheared using various enzyme-based methods. Optimization of the shearing conditions, for example enzyme concentration and incubation time, is needed in order to use enzyme-based methods.

1. DNA End Polishing
 - a. Prepare end repair reaction in a 0.2 ml PCR tube according to Table 1:

Table 1. End Polishing Reaction

Component	Volume
Fragmented DNA	10 μ l (10-50 ng)
10X End Polishing Buffer	1.5 μ l
End Polishing Enzyme Mix	1 μ l
End Polishing Enhancer	1 μ l
Distilled Water	1.5 μ l
Total Volume	15 μl

- b. Mix and incubate for 20 min at 25°C and 20 min at 72°C in a thermocycler (without heated lid).

Note: the amount of fragmented DNA can be 0.2-100 ng with an optimal amount of 10-50 ng.

2. Adaptor Ligation

- a. Prepare a reaction mix for adaptor ligation according to Table 1. Add the following reagents to a 0.2 ml PCR tube containing end polished DNA from step 1.

Table 2. Adaptor Ligation

Component	Volume
End polished DNA (from step 1)	15 μ l
2X Ligation Buffer	17 μ l
T4 DNA Ligase	1 μ l
Adaptors	1 μ l
Total Volume	34 μl

- b. Mix and incubate for 15 min at 25°C in a thermocycler (without heated lid).

Note: (1) The pre-annealed adapters included in the kit are suitable for both non-barcoded (singleplexed) and barcoded (multiplexed) DNA library preparation and are fully compatible with Illumina platforms, such as MiSeq® or HiSeq™ sequencers. (2) If using adaptors from other suppliers (both single-end and barcode adaptors), make sure they are compatible with Illumina platforms and add the correct amount (final concentration 1.5-2 μ M, or according to the supplier's instruction).

3. Size Selection/Clean-up

3.1. Size selection of Ligated DNA

Note: If the starting DNA amount is less than 50 ng, size selection is not recommended and alternatively, clean-up of ligated DNA can be performed prior to PCR amplification according to step 3.2 of the protocol.

- a. Resuspend **MQ Binding Beads** by vortex.
- b. Add 14 μ l of resuspended **MQ Binding Beads** to the tube of ligation reaction. Mix well by pipetting up and down at least 10 times.
- c. Incubate for 5 minutes at room temperature.
- d. Put the tube on an appropriate magnetic stand until the solution is clear (about 4 minutes). Carefully transfer the supernatant containing DNA to a new tube (**Caution: do not discard the supernatant**). Discard the beads that contain the unwanted large fragments.

- e. Add 10 μ l resuspended beads to the supernatant, mix well and incubate for 5 minutes at room temperature.
- f. Put the PCR tube on an appropriate magnetic stand until the solution is clear (about 4 minutes). Carefully remove and discard the supernatant. **Be careful not to disturb or discard the beads that contain DNA.**
- g. Keep the PCR tube in the magnetic stand and add 200 μ l of freshly prepared 80% ethanol to the tube. Incubate at room temperature for 1 min, and then carefully remove and discard the ethanol.
- h. Repeat Step 3.1g one time, for total of two washes.
- i. Open the PCR tube cap and air dry beads for 10 minutes while the tube is on the magnetic stand.
- j. Resuspend the beads in 12 μ l **Elution Buffer**, and incubate at room temperature for 2 minutes to release the DNA from the beads.
- k. Capture the beads by placing the tube in the magnetic stand for 4 minutes or until the solution is completely clear.
- l. Transfer 11 μ l to a new 0.2 ml PCR tube for PCR amplification.

3.2. Clean-up of Ligated DNA (Optional)

- a. Resuspend MQ Binding Beads by vortex.
- b. Add 34 μ l of resuspended beads to the PCR tube of ligation reaction. Mix thoroughly on a vortex mixer or by pipetting up and down at least 10 times.
- c. Incubate for 5 minutes at room temperature to allow DNA to bind to beads.
- d. Put the PCR tube on an appropriate magnetic stand until the solution is clear (about 4 minutes). Carefully remove and discard the supernatant. Be careful not to disturb or discard the beads that contain DNA.
- e. Keep the PCR tube in the magnetic stand and add 200 μ l of freshly prepared 80% ethanol to the tube. Incubate at room temperature for 1 min, and then carefully remove and discard the ethanol.
- f. Repeat Step 3.2e two times for total of three washes.

- g. Open the PCR tube cap and air dry beads for 10 minutes while the tube is on the magnetic stand.
- h. Resuspend the beads in 12 μ l Elution Buffer, and incubate at room temperature for 2 minutes to release the DNA from the beads.
- i. Capture the beads by placing the tube in the magnetic stand for 4 minutes or until the solution is completely clear.
- j. Transfer 11 μ l to a new 0.2 ml PCR tube for PCR amplification.

4. Library Amplification

a. Prepare the PCR Reactions

Thaw all reaction components including master mix, DNA/RNA free water, primer solution and DNA template. Mix well by vortexing briefly. Keep components on ice while in use, and return to -20°C immediately following use. Add components into each PCR tube/well according to the following table:

Component	Size (μ l)
HiFi Master Mix (2X)	12.5 μ l
Primer U	1 μ l
Primer I (or barcode)	1 μ l
Adaptor Ligated DNA	10.5 μ l
Total Volume	25 μl

Important Note: Use of **Primer I** included in the kit will generate a singleplexed library. For multiplexed library preparation, replace **Primer I** with one of the 12 different barcodes (indexes) contained in the AMPINEXT™ NGS Barcode (Index) Set-12 (Cat. No. ENZ-GEN507). You can also add user-defined barcodes (Illumina compatible) instead of **Primer I**.

b. Program the PCR Reactions

Place the reaction plate in the instrument and set the PCR conditions as follows:

Cycle Step	Temp	Time	Cycle
<i>Activation</i>	98°C	30 sec	1
<i>Cycling</i>	98°C 55°C 72°C	20 sec 20 sec 20 sec	Variable*
<i>Final Extension</i>	72°C	2 min	1

* PCR cycles may vary depending on the input DNA amount. In general, use 10 PCR cycles for 100 ng, 12 cycles for 50 ng, 16 cycles for 5 ng, 18 cycles for 1 ng and 22 cycles for 0.2 ng DNA input. Further optimization of PCR cycle number may be required.

5. Clean-up of Amplified Library DNA

- a. Resuspend **MQ Binding Beads** by vortex.
- b. Add 25 µl of resuspended beads to the PCR tube of amplification reaction. Mix thoroughly on a vortex mixer or by pipetting up and down at least 10 times.
- c. Incubate for 5 minutes at room temperature to allow DNA to bind to beads.
- d. Put the PCR tube on an appropriate magnetic stand until the solution is clear (about 4 minutes). Carefully remove and discard the supernatant. **Be careful not to disturb or discard the beads that contain DNA.**
- e. Keep the PCR tube in the magnetic stand and add 200 µl of freshly prepared 80% ethanol to the tube. Incubate at room temperature for 1 min, and then carefully remove and discard the ethanol.
- f. Repeat Step 7e two times for a total of three washes.
- g. Open the PCR tube cap and air dry beads for 10 minutes while the tube is on the magnetic stand.
- h. Resuspend the beads in 22 µl **Elution Buffer**, and incubate at room temperature for 2 minutes to release the DNA from the beads.
- i. Capture the beads by placing the tube in the magnetic stand for 4 minutes or until the solution is completely clear.
- j. Transfer 20 µl to a new 0.2 ml PCR tube.

Quality of the prepared library can be assessed using an Agilent Bioanalyzer or comparable method. Library fragments should have the correct size distribution (ex: 300 bps at peak size) without adaptors or adaptor-dimers. To check the size distribution, dilute library 5-fold with water and apply it to an Agilent high sensitivity chip. If there is presence of <150 bp adaptor dimers or of larger fragments

than expected, they should be removed. To remove fragments below 150 bps use 0.8X **MQ Binding Beads** (ex: add 16 μ l of **MQ Binding Beads** to 20 μ l of sample) according to steps a-i of section **5. Clean-up of Amplified Library DNA**. To remove fragments above 500 bps follow steps a-l of section **3.2 Size Selection of Ligated DNA**.

The prepared DNA library can be quantified with various DNA library quantification methods.

The prepared library DNA can be stored at -20°C until ready to use for sequencing.

TROUBLESHOOTING GUIDE

Problem	Possible Cause	Suggestion
Low yield of library	Insufficient amount of starting DNA.	To obtain the best results, the amount of input DNA should be >10 ng.
	Insufficient purity of starting DNA.	Ensure that RNA is removed by Rnase treatment before starting library preparation protocol.
	Improper reaction conditions at each reaction step.	Check if the reagents are properly added and incubation temperature and time are correct at each reaction step including End Polishing, Adaptor Ligation, size Selection and Amplification.
	Improper storage of the kit	Ensure that the kit has not exceeded the expiration date. Standard shelf life, when stored properly, is 6 months from date of receipt.
Unexpected peak size of Agilent bioanalyzer trace: Presence of <150 bp adaptor dimers or presence of larger fragments than expected.	Improper ratio of MQ Binding Beads to DNA volume in size selection	Check if the correct volume of MQ Binding Beads is added to the DNA solution accordingly. Proper ratios should remove the fragments of unexpected peak size
	Insufficient ligation	Too much and too little input DNA may cause insufficient ligation, which can shift the peak size of the fragment population to be shorter or larger than expected. Make sure that the ligation reaction is properly processed with proper amount of input DNA.
	Over-amplification of library	PCR artifacts from over-amplification of the library may cause the fragment population to shift higher than expected. Make sure to use proper PCR cycles to avoid this problem.



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com