

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

A complete set of reagents required for carrying out a successful ChIP-Seq starting from mammalian cells or tissues.

The AMPINEXT™ ChIP-Seq High-Sensitivity Kit (Illumina) is a complete set of reagents required for carrying out a successful ChIP-Seq starting from mammalian cells or tissues. The kit is designed to selectively enrich a chromatin fraction containing specific DNA sequences from various species, particularly mammals, and to prepare a ChIP-Seq library for next generation sequencing using Illumina platforms such as Illumina Genome Analyzer II, HiSeq and MiSeq systems. The optimized protocol and components of the kit allow capture of low abundance protein/DNA complexes with minimized non-specific background levels and the ability to construct both non-barcoded (singleplexed) and barcoded (multiplexed) ChIP-Seq libraries quickly with reduced bias.

Ordering Information

[Order Online »](#)

ENZ-GEN509-0012	12Reactions
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

- Optimized buffers and protocol allow minimal ChIP background.
- Increased antibody selectivity and capture efficiency.
- Highly efficient enrichment of targeted DNA. Enrichment ratio of positive to negative control > 500.
- High sensitivity and flexibility: Use for both non-barcoded (singleplexed) and barcoded (multiplexed) DNA library preparation
- Fast and streamlined procedure: the procedure from cell/tissues to library DNA is less than 7 hours.
- Highly convenient: the kit contains all required components for each step of ChIP-Seq.
- Minimized bias: Ultra HiFi amplification and optional PCR-free step.

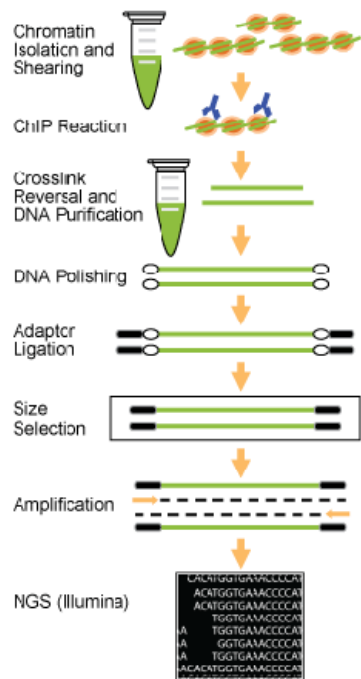


Figure 1. Schematic Procedure for using the AMPINEXT ChIP-Seq High-Sensitivity Kit (Illumina)

Handling & Storage

Use/Stability

Upon receipt: Store the following components at -20°C immediately: Enrichment Enhancer, RNase A, 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: WB, AB, CB, BS, Protease Inhibitor Cocktail, Non-Immune IgG, Anti-RNA Polymerase II, Proteinase K, GAPDH Primer – Forward, GAPDH Primer – Reverse, and 8-Well Assay Strips (With 1 Frame) and 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 reagents required for carrying out a successful ChIP-Seq starting from mammalian cells or tissues

Assay Time

<7 hours

Contents

For ChIP Reaction

WB (Wash Buffer)

AB (Antibody Buffer)

LB (Lysis Buffer)

CB (ChIP Buffer)

DRB (DNA Release Buffer)

DBS (DNA Binding Solution)

BS (Blocker Solution)

DEB (DNA Elution Buffer)

Enrichment Enhancer

Protease Inhibitor Cocktail (PIC)

Non-Immune IgG (1 mg/ml)

Anti-RNA Polymerase II (1 mg/ml)

Proteinase K (10 mg/ml)

RNase A (10 mg/ml)

GAPDH Primer – Forward (20 µM)

GAPDH Primer – Reverse (20 µM)

8-Well Assay Strips (With 1 Frame)

8-Well Strip Caps

Adhesive Covering Film Strip

F-Spin Column

F-Collection Tube

For Library Preparation

10X End Polishing Buffer

End Polishing Enzyme Mix

End Polishing Enhancer

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: March 5, 2025



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