

AMPINEXT™ 5-mC RNA Bisulfite-Seq Easy Kit (Illumina)

A complete set of optimized reagents designed for easily carrying out RNA

The AMPINEXT™ 5-mC RNA Bisulfite-Seq Easy Kit (Illumina) is a complete set of optimized reagents designed for easily carrying out RNA bisulfite conversion, followed by a "post-bisulfite" library preparation process for Illumina platform-based bisulfite sequencing, all in one kit.

Intended applications include whole transcriptome RNA bisulfite sequencing and various other RNA bisulfite-based next generation sequencing techniques for 5-methylcytosine (5-mC) based RNA methylation analysis. The optimized protocol and components of the kit allow the RNA to be bisulfite converted and fragmented simultaneously, followed by quick non-barcoded (singleplexed) and barcoded (multiplexed) library construction using low-nanogram quantities of bisulfite converted RNA.

Manuals, SDS & CofA

[View Online »](#)

- **High sensitivity and efficiency:** Use for non-barcoded (singleplexed) and barcoded (multiplexed) library preparation.
- **Easy and streamlined procedure:** The entire procedure can be finished in 6 hours.
- **Complete conversion:** Converts unmethylated cytosine into uracil at a level greater than 99.9%.
- **Extremely convenient:** The kit contains all the required components for each step of the RNA library preparation process.
- **Minimal bias:** Ultra HiFi amplification.
- **Broad sample suitability:** Starting materials can be total RNA isolated from various tissue/cell samples.

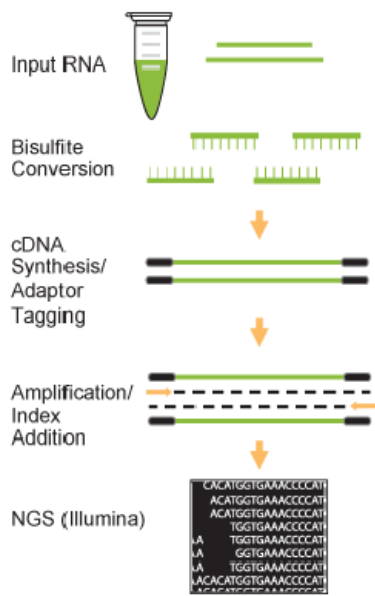


Figure 1. Schematic Procedure for using the AMPINEXT 5-mC RNA Bisulfite-Seq Easy Kit (Illumina)

Handling & Storage

Use/Stability Upon receipt: (1) Store the following components at -20°C immediately: 5X Reaction Buffer, Reaction Enzyme Mix, Adaptor-A, Adaptor-B, RNA Digestion Buffer, RNA Digestion Enzyme, 2X HiFi PCR Master Mix, Primer U, Primer I, and Elution Buffer. (2) Store the following components at 4°C: MQ Binding Beads. (3) Store all other components at room temperature away from light.

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes A complete set of optimized reagents designed for easily carrying out RNA bisulfite conversion, followed by a “post-bisulfite” library preparation process for Illumina platform-based bisulfite sequencing, all in one kit.

Assay Time 6 hours

Contents

- Conversion Buffer
- Conversion Powder
- NA Binding Solution
- F-Spin Column
- F-Collection Tube
- Desulphonation Solution
- 5X Reaction Buffer
- Reaction Enzyme Mix
- Adaptor-A (10 µM)
- Adaptor-B (10 µM)
- RNA Digestion Buffer
- RNA Digestion Enzyme
- MQ Binding Beads
- 2X HiFi PCR Master Mix
- Primer U (10 µM)
- Primer I (10 µM)
- Elution Buffer

Last modified: May 29, 2024



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)