

AMPINEXT™ RNA Bisulfite-Seq Kit (Illumina)

A complete set of optimized reagents designed to carry 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 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.

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------------|
| ENZ-GEN510-0012 | 12Reactions |
| ENZ-GEN510-0024 | 24Reactions |

Manuals, SDS & CofA

[View Online »](#)

- **Fast 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%.
- **High sensitivity, efficiency and flexibility:** Use for both non-barcoded (singleplexed) and barcoded (multiplexed) library preparation.
- **Extremely convenient:** The kit contains all the required components for each step of the RNA library preparation process.
- **Minimal bias:** Ultra HiFi amplification.

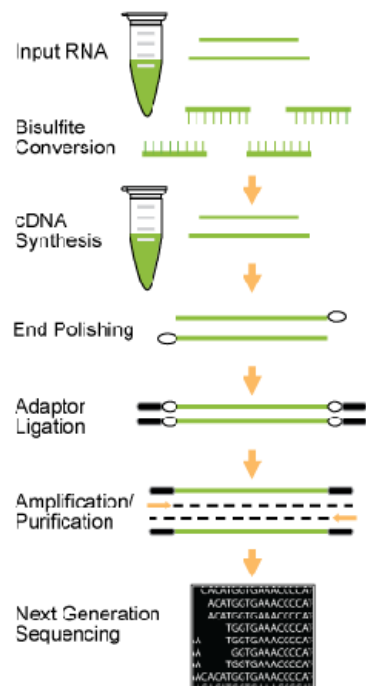


Figure 1. Schematic Procedure for using the AMPINEXT RNA Bisulfite-Seq Kit (Illumina)

Handling & Storage

| | |
|---------------|--|
| Use/Stability | Upon receipt: (1) Store the following components at -20°C immediately: 5X RT Reaction Buffer, 10 mM dNTP mix, 0.1M DTT, RNase Inhibitor, Random Primer, RT Enzyme Mix, 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. |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|-------------------|---|
| Application Notes | A complete set of optimized reagents designed to carry 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 RT Reaction Buffer 10 mM dNTP Mix 0.1M DTT RNase Inhibitor Random Primer (50 µM) RT Enzyme Mix 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 |



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)