

# 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."

The AMPINEXT™ RNA Bisulfite-Seq Kit (Illumina) is 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
-----------------	-------------

## 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.



## 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

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)