

Instruction Manual

Ribominus™ Transcriptome Isolation Kit (Yeast and Bacteria)

For efficient transcriptome enrichment by depleting large ribosomal RNA from yeast and bacteria total RNA

Catalog nos. K1550-03, K1550-04, K1550-05

Version A

12 September 2005
25-0879

Table of Contents

Table of ContentsIII
Experienced Users Procedurev
Kit Contents and Storagevii
Accessory Productsx
Overview1
Product Specifications4
Experimental Overview6
Preparing Total RNA7
Selective Hybridization and Removal of rRNA9
Using RiboMinus [™] Concentration Module to Concentrate RiboMinus [™] RNA17
Using Ethanol Precipitation to Concentrate RiboMinus [™] RNA20
Analyzing RiboMinus [™] RNA21
Troubleshooting24
Technical Service26
Purchaser Notification27
References28

Experienced Users Procedure

Introduction

This quick reference protocol is included for experienced users of the RiboMinusTM Yeast Transcriptome Isolation Kit and RiboMinusTM Bacteria Transcriptome Isolation Kit. If you are a first time user, follow the detailed protocol in this manual.

Step	Action			
Isolating Total RNA	Isolate high-quality total RNA from yeast or bacteria cells using a method of choice prior to using these kits.			
	You will need 2-10 μg total RNA in <20 μl per reaction.			
Preparing RiboMinus™	Resuspend the RiboMinus [™] Magnetic Beads in its bottle by thorough vortexing.			
Magnetic Beads	Pipet 250 μ l of the bead suspension into a sterile, RNase-free, 1.5-ml microcentrifuge tube.			
	3. Place the tube with the bead suspension on a magnetic stand for 1 minute. Gently aspirate and discard the supernatant.			
	4. Add 250 μl sterile, RNase-Free Water to the beads and resuspend beads. Place the tube on a magnetic stand for 1 minute. Gently aspirate and discard the supernatant.			
	5. Repeat Step 4 once.			
	6. Resuspend beads in 250 μ l Hybridization Buffer (B10). Place the tube on a magnetic stand for 1 minute. Gently aspirate and discard the supernatant.			
	7. Resuspend beads in 100 μl Hybridization Buffer (B10) and keep the beads at room temperature until use.			
Selective Hybridization	Perform hybridization of your total RNA sample with the RiboMinus [™] Yeast Probe or RiboMinus [™] Bacteria Probe, depending on the source of total RNA as below.			
	1. To a sterile, RNase-free 1.5 ml microcentrifuge tube, add:			
	Total RNA (2-10 μg): <20 μl			
	RiboMinus™ Probe (100 pmol/μl): 4 μl			
	Hybridization Buffer (B10): 100 μl			
	2. Incubate the tube at 37°C for 5 minutes to denature RNA.			
	3. Place the sample on ice for at least 30 seconds			
	4. Proceed to Removing rRNA, next page.			

Experienced Users Procedure, Continued

Step	Action
Removing rRNA	Briefly centrifuge the tube with the cooled hybridized sample (from Step 3, previous page) to collect the sample to the bottom of the tube.
	 Transfer ~124 μl of the cooled hybridized sample (from Step 3, previous page) to the prepared RiboMinus[™] Magnetic beads from Step 7, previous page, and mix well by vortexing.
	3. Incubate the tube at 37°C for 15 minutes. During incubation, gently mix the contents occasionally.
	4. Place the tube on a magnetic stand for 1 minute to pellet the rRNA-probe complex. The supernatant contains the RiboMinus™ RNA fraction.
	5. Transfer the supernatant (~224 μl) to a tube capable of holding 3X volume of the supernatant.
Concentrating RiboMinus [™] RNA using	See below for concentrating the RiboMinus [™] RNA using the RiboMinus [™] Concentration Module (pages 17-19) or using ethanol precipitation (page 20).
RiboMinus [™] Concentration Module	1. To the sample from Step 5, above, add 250 μl Binding Buffer (L3) and 125 μl 96-100% ethanol. Mix well by vortexing.
	2. Bind the sample from Step 1 containing Binding Buffer (L3) and ethanol to the spin column. Centrifuge the column at $12,000 \times g$ for 1 minute at room temperature. Discard the flow through.
	3. Wash the column with 200 µl Wash Buffer (W5) with ethanol (page 18). Centrifuge the column at ≥12,000 × g for 1 minute at room temperature. Discard the flow through.
	4. Repeat the wash step once.
	5. Discard the tube and place the column into a clean Wash Tube, supplied with the kit.
	6. Centrifuge the column at ≥12,000 × g for 2–3 minutes at room temperature to remove any residual Wash Buffer (W5). Place the column in a 1.5-ml Recovery Tube.
	7. Elute with 10-15 μl of RNase-Free Water. Incubate the column at room temperature for 1 minute. Centrifuge the column at ≥12,000 × g for 1 minute.
	The Recovery Tube contains purified RiboMinus™ RNA.
	8. Store RiboMinus™ RNA at -80°C or place RiboMinus™ RNA on ice to proceed to desired downstream application.

Kit Contents and Storage

Types of Kits

This manual is supplied with the following products.

Product	Catalog no.
RiboMinus [™] Yeast Transcriptome	K1550-03
Isolation Kit	
RiboMinus™ Bacteria Transcriptome	K1550-04
Isolation Kit	
RiboMinus [™] Concentration Module	K1550-05

Shipping and Storage

All components of the RiboMinus $^{\text{\tiny TM}}$ Transcriptome Isolation Kits and the RiboMinus $^{\text{\tiny TM}}$ Concentration Module are shipped at room temperature.

Upon receipt, store as follows:

- RiboMinus[™] Transcriptome Isolation Kits at 4°C
- RiboMinus[™] Concentration Module at room temperature

RiboMinus[™] Transcriptome Isolation Kits

The components included in the RiboMinus $^{\text{\tiny TM}}$ Transcriptome Isolation Kits are listed below.

Sufficient reagents are provided in the kits to perform 12 reactions.

Store the module at 4°C. For long-term storage, store the probe at -20°C.

Note: Some reagents are provided in excess.

Component	K1550-03	K1550-04
RiboMinus [™] Magnetic Beads (12 mg/ml) in Phosphate Buffered Saline (PBS), pH 7.4 containing 0.01% Tween 20 and 0.09% sodium azide	3 ml	3 ml
RiboMinus [™] Yeast Probes in ultrapure water (100 pmol/µl)	50 μl	-
RiboMinus™ Bacteria Probes in ultrapure water (100 pmol/µl)	-	50 μ1
Hybridization Buffer (B10)	6 ml	6 ml
RNase-Free Water	6 ml	6 ml

Kit Contents and Storage

RiboMinus[™] Concentration Module

The components included in the RiboMinus $^{\text{\tiny M}}$ Concentration Module are listed below.

All components are shipped at room temperature.

Sufficient reagents are provided in the kit to perform 6 reactions.

Store the Concentration Module at room temperature.

Note: Some reagents are provided in excess.

Component	Amount
Binding Buffer (L3)	3.3 ml
Wash Buffer (W5)	1.5 ml
RNase-Free Water	6.6 ml
Spin Columns with Collection Tubes	6
Wash Tubes (2.0 ml)	6
Recovery Tubes (1.5 ml)	6



The reagents provided in the RiboMinus[™] Yeast or Bacteria Transcriptome Isolation Kits are sufficient for twice as many reactions (12 reactions) as provided with the RiboMinus[™] Concentration Module (6 reactions).

To concentrate RiboMinus[™] RNA purified with RiboMinus[™] Yeast or Bacteria Transcriptome Isolation Kits with the RiboMinus[™] Concentration Module, order **two** RiboMinus[™] Concentration Modules per RiboMinus[™] Yeast or Bacteria Transcriptome Isolation Kit.

Kit Contents and Storage, Continued

Product Qualification

The RiboMinus[™] Yeast Transcriptome Isolation Kit and RiboMinus[™] Bacteria Transcriptome Isolation Kit are functionally qualified as described below.

Purified total RNA (10 μ g) from yeast (*S. cerviseiae*) is subjected to ribosomal RNA depletion using the RiboMinus[™] Yeast Transcriptome Isolation Kit as described in this manual.

Purified total RNA (10 μ g) from bacteria *E. coli* (TOP10) is subjected to ribosomal RNA depletion using the RiboMinus[™] Bacteria Transcriptome Isolation Kit as described in this manual.

Agarose gel electrophoresis must show >98% depletion of the large ribosomal RNA bands (yeast: 18S and 25/26S rRNA; bacteria: 16S and 23S rRNA) from the purified sample as compared to control samples.

Bioanalyzer analysis of the purified sample must show <2% of peak area for the large ribosomal bands as compared to peak area for the control sample.

RiboMinus™ Magnetic Beads

The binding capacity of the beads must be >2500 pmoles free biotin per mg of streptavidin-coated magnetic beads and must be free from bacterial contamination.

RiboMinus[™] Probes

The probes must contain the correct sequence and the locked nucleic acid (LNA $^{\text{\tiny TM}}$) at the specified position for each probe. Mass spectrometry analysis of probes must indicate the specified mass and HPLC analysis must indicate >85% purity. The probes must be RNase- and DNase-free.

Purchaser Notification

Limited Use Label License No: 237 LNA™ Oligonucleotides

 $LNA^{\text{\tiny TM}}$ oligonucleotides are produced under a license from Exigon A/S.

Accessory Products

Additional Products

The following products are also available from Invitrogen. For more details on these products, visit our Web site at www.invitrogen.com or contact Technical Service (page 26).

Product	Quantity	Catalog no.
RNase AWAY®	250 ml	10328-011
UltraPure™ DEPC-treated Water	1 L	750023
UltraPure™ DNase/RNase-Free Distilled Water	500 ml	10977-015
Quant-iT™ RNA Assay Kit	1000 assays	Q-33140
Micro-to-Midi™ Total RNA Purification System	50 reactions	12183-018
TRIzol® Reagent	100 ml	15596-026
DNase I	20,000 units	18047-019
DNase I, Amplification Grade	100 units	18068-015
Magna-Sep [™] Magnetic Particle Separator	1	K1585-01
RiboMinus™ Human/Mouse Transcriptome Isolation Kit	1	K1550-02

Overview

Introduction

The RiboMinus[™] Transcriptome Isolation Kits provide a novel and efficient method to isolate RNA molecules of the transcriptome devoid of large ribososmal RNA (rRNA) from total RNA for transcriptome analysis. The purification method is not dependent on the polyadenylation status or presence of a 5′-cap structure on the RNA. See below for more details on the purification protocol.

The isolation of RNA fraction depleted of ribososmal RNA is achieved by the selective removal of the large rRNA molecules from total RNA. The resulting rRNA depleted RNA fraction is termed as RiboMinus™ RNA (see next page for details).

Using the kit to isolate RiboMinus™ RNA results in efficient (>98%) removal of large rRNA molecules (yeast: 18S and 25/26S; bacteria 16S and 23S) from 10 µg total RNA enabling the analysis of the whole transcriptome without any interference from rRNA that account for ~90-95% RNA species in total RNA.

System Overview

The RiboMinus™ Transcriptome Isolation Kits are based on the selective removal of abundant large ribosomal RNA molecules from total RNA and concentrating the RiboMinus™ RNA enriched fraction.

Total RNA is hybridized with species-specific, rRNA sequence-specific 5′-biotin labeled oligonucleotide probes (RiboMinus[™] Yeast Probes or RiboMinus[™] Bacteria Probes). The rRNA/5′-biotin labeled probe complex is removed from the sample with streptavidin coated magnetic beads (RiboMinus[™] Magnetic Beads).

The RiboMinus[™] RNA sample is then concentrated using RiboMinus[™] Concentration Module for a spin column-based centrifugation protocol (page 17) or using ethanol precipitation (page 20). The binding conditions of the spin column method are optimized for the RiboMinus[™] RNA sample with ethanol and Binding Buffer (L3). The sample is loaded onto a spin column. The RiboMinus[™] RNA binds to the silica-based membrane in the column and impurities are removed by thorough washing with Wash Buffer (W5). The RNA is then eluted in sterile RNase free water.

For details on RiboMinus[™] Probes and RiboMinus[™] Magnetic Beads, see page 9.

Overview, Continued

Advantages

Using the RiboMinus[™] Transcriptome Isolation Kits to isolate RiboMinus[™] RNA (rRNA depleted RNA) provides the following advantages:

- Rapid and efficient isolation of high-quality RiboMinus[™] RNA using probes specific to the large rRNA species
- Specifically designed to isolate RiboMinus[™] RNA enriched in (polyA) mRNA (yeast), nonpolyadenylated RNA, pre-processed RNA, tRNA, and small rRNAs
- Minimal contamination from rRNA molecules
- Reliable performance of the RiboMinus[™] RNA in downstream applications such as microarray analysis, cDNA library construction, and qRT-PCR

RiboMinus[™] RNA

The large ribososmal RNA depleted RNA fraction is termed as RiboMinus^T RNA fraction.

The RiboMinus[™] RNA fraction contains polyadenylated (polyA) mRNA (yeast), non-polyadenylated RNA, pre-processed RNA, tRNA, small rRNAs (5S rRNA, and additional 5.8S rRNA for eucaryotic RNA), and may also contain regulatory RNA molecules such as microRNA (miRNA) and short interfering RNA (siRNA), snRNA, and other RNA transcripts of yet unknown function.

The RiboMinus™ RNA molecules are part of the transcriptome and are important in protein coding, signaling, structural support of subcellular elements, and transcriptional/post transcriptional regulation.

The transcriptome is defined as the complete collection of transcribed elements of the genome (Ruan *et al.*, 2004) and contains mRNA transcripts and non-mRNA transcripts including RiboMinus[™] RNA. Transcriptome analysis is gaining increased attention in gene expression analysis. Since large rRNA constitutes 90-95% RNA species in total RNA, whole transcriptome analysis without any contamination from rRNA is very difficult and suggests the need for developing procedures for transcriptome isolation.

Overview, Continued

Drawbacks of RNA Purification Methods

Current methods for RNA purification do not allow for efficient isolation of transcriptome. The total RNA purification methods result in enriching the large rRNA molecules while the mRNA purification methods use polyA-selection and/or cap-binding approaches (for eucaryotic RNA) that do not enrich the complete transcriptome.

The RiboMinus[™] Transcriptome Isolation Kits offer a novel method of isolating transcriptome and involves selective removal of large rRNA from total RNA. The isolated transcriptome is >98% depleted in rRNA and is enriched in all RNA transcripts of interest enabling whole transcriptome analysis.

Downstream Applications

The isolated RiboMinus[™] RNA is suitable for use in downstream applications such as microarray analysis, qRT-PCR, and cDNA library construction

Product Specifications

System Specifications Starting Material: 2-10 μg total RNA in <20 μl

rRNA Removal: >98%

RiboMinus[™] RNA Yield: ≥0.5 µg from 10 µg total RNA

RiboMinus[™] Probes Specifications RiboMinus[™] Yeast Probes:

Probe Contents: 2 probes each for 18S and

25/26S rRNA

Probe Specificity: Yeast

Probe Size: 18-21 oligonucleotides

Probe Label: 5'-biotin label

LNA[™] Content: Each probe contains 5-7

LNA[™] monomers in the

oligonucleotide

Probe Mixture Concentration: 100 pmol/μl

RiboMinus[™] Bacteria Probes:

Probe Contents: 2 probes each for 16S and

23S rRNA

Probe Specificity: Bacteria

Probe Size: 19-20 oligonucleotides

Probe Label: 5'-biotin label

LNA[™] Content: Each probe contains 4-5

LNA[™] monomers in the

oligonucleotide

Probe Mixture Concentration: 100 pmol/µl

For details on the probe, see page 9.

RiboMinus[™] Magnetic Bead Specifications The RiboMinus $^{\scriptscriptstyle{\text{TM}}}$ Magnetic Beads are streptavidin-coated

magnetic beads.

Bead Binding Capacity: >2500 pmoles free biotin per mg

RiboMinus[™] Magnetic Beads

Bead Size: 1 µm diameter

Magnet Particle: Superparamagnetic polydisperse

core-shell polystyrene particles

Concentration: 12 mg/mlSpecific Gravity: $1.1-1.4 \text{ g/cm}^3$ For details on the beads, see page 10.

Product Specifications, Continued

RiboMinus[™] Concentration Module Specifications

Binding Capacity: ~5 μg nucleic acid

Column Reservoir Capacity: $700 \ \mu l$ Wash Tube Capacity: $2.0 \ ml$ Recovery Tube Capacity: $1.5 \ ml$

Centrifuge Compatibility: Capable of centrifuging at

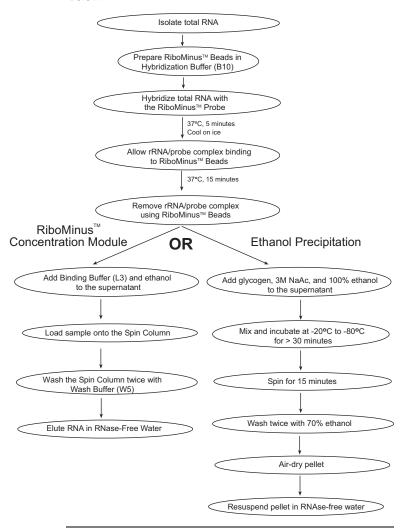
 $>10,000 \times g$

Elution Volume: $10-15 \mu l$

Experimental Overview

Introduction

The flow chart for isolating transcriptome using the RiboMinus $^{\text{\tiny TM}}$ Yeast Transcriptome Isolation Kit and RiboMinus $^{\text{\tiny TM}}$ Bacteria Transcriptome Isolation Kit is shown below.



Preparing Total RNA

Introduction

You will need to isolate high-quality total RNA from cells or tissues using a method of choice prior to using this kit.

To obtain high-quality total RNA, follow the guidelines recommended below.

General Handling of RNA

Observe the following guidelines to prevent RNase contamination:

- Use disposable, individually wrapped, sterile plasticware
- Use only sterile, new pipette tips and microcentrifuge tubes
- Wear latex gloves while handling reagents and RNA samples to prevent RNase contamination from the surface of the skin
- Always use proper microbiological aseptic techniques when working with RNA
- Use RNase AWAY® Reagent (page x) to remove RNase contamination from surfaces

Total RNA Isolation

Total RNA can be isolated from tissue or cells using the method of choice. We recommend isolating total RNA using the Micro-to-Midi $^{\text{\tiny{M}}}$ Total RNA Purification System or TRIzol $^{\text{\tiny{B}}}$ Reagent available from Invitrogen (see page x for ordering information).

You will use 2-10 μ g total RNA in less than 20 μ l for each reaction. Resuspend isolated total RNA in DEPC-treated water accordingly (\geq 0.5 μ g/ μ l).

Check the quality of your total RNA, including DNA contamination (see below). Store your total RNA at -80°C and avoid repeated freezing and thawing of total RNA.



If your downstream application requires DNA-free RNA, perform DNase-treatment of the total RNA **before** purifying RiboMinus $^{\text{\tiny{TN}}}$ RNA.

Preparing Total RNA, Continued

Checking the Total RNA Quality

To check total RNA integrity, analyze $\sim\!0.5~\mu g$ of your RNA by agarose/ethidium bromide gel electrophoresis. You should see the following on an agarose gel:

- for yeast total RNA: 18S rRNA band (~2.0 kb) and 25/26S rRNA band (~3.8 kb)
- for bacterial total RNA: 16S rRNA band (~1.5 kb) and 23S rRNA band (~2.9 kb)
- The larger bands should be approximately twice the intensity of the smaller bands

Selective Hybridization and Removal of rRNA

Introduction

Instructions are provided in this section for selective hybridization of rRNA to the RiboMinus[™] Probe and removal of rRNA using RiboMinus[™] Magnetic Beads. See page 11 for an experimental outline.

RiboMinus[™] Probes

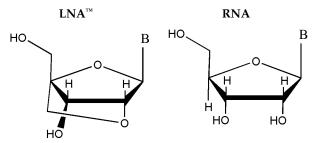
The RiboMinus™ Yeast Probes and RiboMinus™ Bacteria Probes are mixtures of oligonucleotide probes containing 2 probes each specific for 18S rRNA and 25/26S rRNA (yeast) or 16S rRNA and 23S rRNA (bacteria), respectively (see page 4 for specifications). The probes are designed to hybridize species-specific with highly conserved regions of the large rRNA molecules.

Each probe is single-stranded and contains 4-7 LNA $^{\text{IM}}$ (\underline{L} ocked \underline{N} ucleic \underline{A} cid) monomers incorporated at specific locations. The incorporation of LNA $^{\text{IM}}$ (see next page for details on LNA $^{\text{IM}}$) into the oligonucleotide probe increases the depletion efficiency of the rRNA from the samples without increasing the amount of beads or probe concentration.

The 5'-end of each probe is conjugated to biotin to allow removal of rRNA/probe complexes by binding to stretptavidin RiboMinus[™] Magnetic Beads (see next page).

LNA[™] (Locked Nucleic Acid)

The structure of the LNA $^{\text{\tiny TM}}(\underline{L}\text{ocked }\underline{N}\text{ucleic }\underline{A}\text{cid})$ monomer (see figure below) consists of a ribonucleoside linked between the 2' oxygen and 4' carbon atom of the methylene ring (Braasch and Corey, 2001).



This configuration locks the sugar backbone resulting in an increase in T_m (melting temperature).

Incorporation of 4-7 LNA $^{\text{\tiny IM}}$ monomers into an oligonucleotide does not affect the ability of the oligonucleotide to bind DNA or RNA but increases the stability of the oligonucleotide/RNA complex (McTigue et al., 2004). Oligonucleotides containing LNA $^{\text{\tiny IM}}$ are used in hybridization assays requiring high specificity and reproducibility.

RiboMinus[™] Magnetic Beads

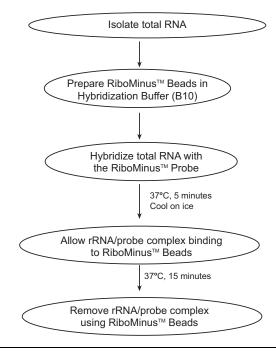
The RiboMinus $^{\text{\tiny{TM}}}$ Magnetic Beads are streptavidin-coated magnetic beads used for the removal of probe/rRNA complexes from the sample. The beads bind to the biotin-labeled probe complexed with rRNA or the probe alone.

The beads are 1 μ m polystyrene beads with a magnetic core that is strong enough to separate the bound complex from the solvent in a short period of time (see page 4 for specifications). The beads do not promote non-specific binding of any other RNA molecules.

The size and the biotin binding capacity of the RiboMinus $^{\text{\tiny M}}$ Magnetic Beads is optimized for use with this kit and results in >98% depletion of rRNA using 10 μg total RNA as the starting material. Avoid using any other streptavidin-coated magnetic beads with this kit.

Experimental Outline

The figure below depicts the experimental outline for hybridization of rRNA to specific probes and removal of rRNA.



Materials Needed

You will need the following items:

- Total RNA (see pages 7-8)
- Magna-Sep[™] Magnetic Particle Separator (page x) or equivalent
- RNase-free microcentrifuge tubes
- Water baths or heat blocks set to 70-75°C and 37°C
- Ice

Components supplied with the kit

- RiboMinus[™] Magnetic Beads, keep on ice until use
- RiboMinus[™] Yeast Probe or RiboMinus[™] Bacteria Probe, keep on ice until use
- Hybridization Buffer (B10)
- RNase-Free Water



Follow the recommendations for handling the RiboMinus™ Magnetic Beads below for best results:

- During the mixing and washing steps of the magnetic beads, mix beads using a vortex. A low speed centrifuge pulse may be required to remove beads stuck in the tube cap. Avoid mixing by pipetting up and down as it results in bead loss.
- Do not allow the beads to dry as drying reduces the bead efficiency. During all washing steps with beads, add water or buffer to the tube containing beads while the tube is still on a magnetic stand to prevent drying of beads. Remove the tube from the magnet and resuspend the beads as described above.
- To aspirate the supernatant after bead washing, place the pipette tip at the opposite side of the tube, away from the beads. Carefully remove the supernatant without disturbing or removing any beads.
- Do not submerge the magnetic stand in water. To clean the magnetic stand, spray the stand with ethanol and wipe it with a paper towel.



The RiboMinus[™] Hybridization Buffer (B10) contains guanidine isothiocyanate. Always wear a laboratory coat, disposable gloves, and eye protection when handling buffers.

Do not add bleach or acidic solutions directly to solutions containing guanidine isothiocyanate or sample preparation waste as it forms reactive compounds and toxic gases when mixed with bleach or acids.

Preparing RiboMinus[™] Magnetic Beads

Follow the recommendations on page 12 for handling beads and performing the washing steps.

- Resuspend the RiboMinus[™] Magnetic Beads in its bottle by thoroughly vortexing.
- 2. Pipet 250 µl of the bead suspension into a sterile, RNase-free, 1.5 ml microcentrifuge tube.
- 3. Place the tube with the bead suspension on a magnetic separator for 1 minute. The beads will settle to the side of the tube that faces the magnet. Aspirate and discard the supernatant.
- 4. Add 250 μl sterile, RNase-Free Water supplied with the kit to the beads and resuspend beads by vortexing.
- 5. Place the tube on a magnetic separator for 1 minute. Gently aspirate and discard the supernatant.
- 6. Repeat Steps 4-5 once.
- Resuspend beads in 250 μl Hybridization Buffer (B10).
 Place the tube on a magnetic separator for 1 minute.
 Gently aspirate and discard the supernatant.
- 8. Resuspend beads in $100 \mu l$ Hybridization Buffer (B10) and keep the beads at 37° C until use

Continue to Hybridization Step, next page.

Hybridization Step

Instructions are provided below to perform hybridization for 2-10 μg of your total RNA sample with the RiboMinusTM Yeast Probe or RiboMinusTM Yeast Probe, depending o the species of the total RNA.

If you wish to process >10 μg total RNA sample, divide your sample into two samples, each containing <10 μg total RNA.

- 1. Set a water bath or heat block to 37°C.
- 2. To a sterile, RNase-free 1.5 ml microcentrifuge tube, add the following:

Total RNA (2-10 μg):	<20 μl
RiboMinus [™] Probe (100 pmol/μl):	4 μl
Hybridization Buffer (B10):	100 μl

- 3. Incubate the tube at 37°C for 5 minutes to denature RNA.
- 4. Place the sample on ice for at least 30 seconds.
- 5. Proceed to Removing rRNA, next page

Removing rRNA

- 1. Set a water bath or heat block to 37°C.
- Briefly centrifuge the tube with the cooled hybridized sample (from Step 4, previous page) to collect the sample to the bottom of the tube.
- Transfer the sample (~124 µl) to the prepared RiboMinus[™]
 Magnetic beads from Step 8, page 14. Mix well by
 vortexing the tube repeatedly.
- 4. Incubate the tube at 37°C for 15 minutes. During incubation, gently mix the contents occasionally. Briefly centrifuge the tube to collect the sample to the bottom of the tube.
- Place the tube on a magnetic separator for 1 minute to pellet the rRNA-probe complex. Do not discard the supernatant. The supernatant contains RiboMinus™ RNA.
- 6. Transfer the supernatant (\sim 224 μ l) to a tube capable of holding 3X the volume of the supernatant.

Next Step

You can concentrate the RiboMinus[™] RNA using RiboMinus[™] Concentration Module with a rapid and optimized spin-based protocol (pages 17-19) or using ethanol precipitation (page 20).

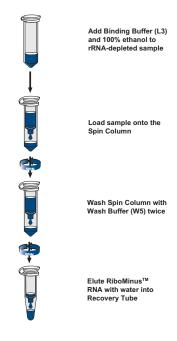
Using RiboMinus[™] Concentration Module to Concentrate RiboMinus[™] RNA

Introduction

The RiboMinus™ Concentration Module is designed for concentrating RiboMinus™ RNA using spin column-based centrifugation protocol in a total time of **10-15 minutes**. See next page for an experimental outline.

Experimental Outline

The figure below depicts the experimental outline for concentrating the RiboMinus $^{\text{\tiny M}}$ RNA using a spin-column based centrifugation procedure.



Using RiboMinus[™] Concentration Module to Concentrate RiboMinus[™] RNA, Continued

Materials Needed

- RiboMinus[™] RNA sample from Step 6, page 16
- 96-100% ethanol
- Microcentrifuge capable of centrifuging >12,000 x g
 Components supplied with the RiboMinus[™] Concentration Module
- Binding Buffer (L3)
- Wash Buffer (W5)
- RNase-Free Water
- Spin Column with Collection Tubes
- Wash Tubes
- Recovery Tubes



The RiboMinus[™] Concentration Module Binding Buffer (L3) contains guanidine isothiocyanate. Always wear a laboratory coat, disposable gloves, and eye protection when handling buffers.

Do not add bleach or acidic solutions directly to solutions containing guanidine isothiocyanate or sample preparation waste as it forms reactive compounds and toxic gases when mixed with bleach or acids.

Before Starting

Add 6 ml 96-100% ethanol to 1.5 ml Wash Buffer (W5) included with the kit. Store the Wash Buffer (W5) with ethanol at room temperature.

Binding Step

- 1. To the sample from Step 6, page 16, add 250 μ l Binding Buffer (L3) and 125 μ l 96-100% ethanol. Mix well by vortexing.
- 2. Load the sample containing Binding Buffer (L3) and ethanol to the column.
- 3. Centrifuge the column at ≥12,000 × g for 1 minute at room temperature. Discard the flow through.
- 4. Proceed to **Washing Step**, next page.

Using RiboMinus[™] Concentration Module to Concentrate RiboMinus[™] RNA, Continued

Washing Step

- 1. Add 200 μl Wash Buffer (W5) with ethanol (page 18) to the column.
- 2. Centrifuge the column at ≥12,000 × g for 1 minute at room temperature. Discard the flow through.
- 3. Repeat the wash step with 200 µl Wash Buffer (W5) with ethanol.
- 4. Discard the collection tube and place the column into a clean Wash Tube supplied with the kit.
- Centrifuge the column at maximum speed for 2-3 minutes at room temperature to remove any residual Wash Buffer (W5). Discard the Wash Tube.
- 6. Proceed to **Elution Step**, below.

Elution Step

- 1. Place the Spin Column in a clean 1.5-ml Recovery Tube supplied with the kit.
- 2. Add 10-15 μ l of RNase-Free Water to the center of the column. Incubate the column at room temperature for 1 minute.
- 3. Centrifuge the column at maximum speed for 1 minute at room temperature.
 - The Recovery tube contains purified and concentrated RiboMinusTM RNA sample that is depleted of rRNA.
- Place RiboMinus[™] RNA on ice to proceed to desired downstream application or store RiboMinus[™] RNA at -80°C until further use.
 - See pages 21-23 to analyze yield and quality of the RiboMinus™ RNA.

Using Ethanol Precipitation to Concentrate RiboMinus[™] RNA

Introduction

This section includes a protocol for **Ethanol Precipitation** to further concentrate the RiboMinus^m RNA. You will need at least 1 hour to perform the ethanol precipitation.

Materials Needed

- RiboMinus[™] RNA sample (Step 7, page 16)
- Glycogen, 20 μg/μl (see page x)
- 3 M sodium acetate in RNAse-free water
- 96-100% cold ethanol
- 70% cold ethanol
- RNase-free water
- Sterile, RNase-free microcentrifuge tubes
- Microcentrifuge capable of centrifuging >12,000 x g

Ethanol Precipitation

- 1. Add the following components to the RiboMinus[™] RNA:
 - 1 μl glycogen (20 μg/μl)
 - 1/10th sample volume of 3 M sodium acetate
 - 2.5X sample volumes of 100% ethanol
- 2. Mix well and incubate at -20 or -80°C for a minimum of 30 minutes.
- 3. Centrifuge the tube for 15 minutes \geq 12,000 × g at 4°C.
- 4. Carefully discard the supernatant without disturbing the pellet.
- 5. Add 500 µl 70% cold ethanol.
- 6. Centrifuge the tube for 5 minutes at \geq 12,000 × g at 4°C.
- 7. Carefully discard the supernatant without disturbing the pellet.
- 8. Repeat Steps 6-8 once.
- 9. Air-dry the pellet for ~5 minutes; do not dry completely
- 10. Resuspend the RNA pellet in ~10-30 μl RNase-free water.
- Place RiboMinus[™] RNA on ice to proceed to desired downstream application or store RiboMinus[™] RNA at -80°C until further use.

See pages 21-23 to analyze yield and quality of RiboMinus™ RNA.

Analyzing RiboMinus[™] RNA

RNA Yield

The quantity of the purified RiboMinus^m RNA is easily quantitated using UV absorbance at 260 nm or Quant-iT^m RNA Assay Kit.

UV Absorbance

1. Dilute an aliquot of the small sample in 10 mM Tris-HCl, pH 7.0. Mix well. Transfer to a cuvette (1-cm path length).

Note: The RNA must be in a neutral pH buffer to accurately measure the UV absorbance.

 Determine the OD₂₆₀ of the solution using a spectrophotometer blanked against 10 mM Tris-HCl, pH 7.0.

Calculate the amount of total RNA using the following formula:

Total RNA (μ g) = OD₂₆₀ x 40 μ g/(1 OD₂₆₀ x 1 ml) x dilution factor x total sample volume (ml)

The typical yield of RNA using the RiboMinus $^{\scriptscriptstyle{\text{TM}}}$ Transcriptome Isolation Kits (Yeast and Bacteria) is $\sim\!0.5~\mu g$ RNA from 10 μg total RNA sample.

Quant-iT™ RNA Assay Kits

The Quant-iT™ RNA Assay Kit (see page x) provides a rapid, sensitive, and specific method for RNA quantitation with minimal interference from DNA, protein, or other common contaminants that affect UV absorbance readings.

The kit contains a state-of-the-art quantitation reagent and pre-diluted standards for standard curve. The assay is performed in a microtiter plate format and is designed for reading in standard fluorescent microplate readers.

Analyzing RiboMinus[™] RNA, Continued

RNA Quality

The RNA isolated using the RiboMinus™ Transcriptome Isolation Kit is of high-quality and is >98% depleted in rRNA species.

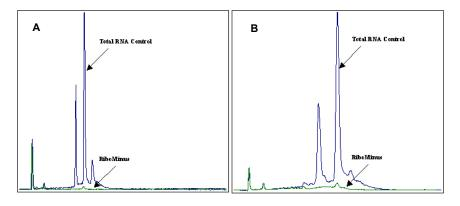
To verify rRNA depletion, perform agarose gel electrophoresis of the sample or use a bioanalyzer (see next page for examples).

Gel Electrophoresis to assess RNA Quality

You can use agarose gel electrophoresis analysis to verifyl depletion of the large rRNA bands as compared to a control sample and absence of contaminating DNA and RNA degradation.

Analyzing RiboMinus[™] RNA, Continued

Bioanalyzer to assess RiboMinus[™] RNA Quality The efficiency of rRNA depletion in RiboMinus[™] RNA, RNA degradation, and RNA concentration can be analyzed using a bioanalyzer such as the Agilent 2100 bioanalyzer with an RNA LabChip[®]. In the example below, the bioanalyzer data was used to show efficient removal of 18S and 25/26S rRNA from yeast total RNA and of 16S and 23S rRNA from bacterial total RNA.



RiboMinus[™] RNA was purified from 10 µg total RNA from *S. cerviseiae* yeast cells (panel A) using RiboMinus[™] Transcriptome Isolation Kit (Yeast) or from *E. coli* bacteria (panel B) using RiboMinus[™] Transcriptome Isolation Kit (Bacteria) as described in this manual. The total RNA control samples were obtained by omitting RiboMinus[™] Magnetic Beads in the reaction mixture. Aliquots of 2% of the final elution volume of RiboMinus[™] RNA and control total RNA were subjected to the bioanalysis using Agilent 2100 bioanalyzer. The graph shows the removal of the large rRNA from the purified RiboMinus[™] RNA samples as compared to control samples.

23

Troubleshooting

Introduction

Review the table below to troubleshoot problems that you may encounter using the RiboMinus $^{\text{\tiny M}}$ Transcriptome Isolation Kits.

Problem	Cause	Solution	
Low RNA yield	Low RNA content	Various tissues have different RNA content and the yield is dependent on the sample.	
	Incorrect binding conditions when using the RiboMinus™ Concentration Module	For efficient binding of RiboMinus [™] RNA to the spin column, always add 250 µl Binding Buffer (L3) and 125 µl 96-100% ethanol to the hybridized sample and mix well by vortexing prior to loading onto the Spin Column.	
	Ethanol not added to Wash Buffer (W5)	Be sure to add 96–100% ethanol to Wash Buffer (W5) as described on page 18.	
	Incorrect elution conditions	Add water to the center of the column and perform incubation for 1 minute with water before centrifugation.	
	RNA quantitation performed with water	Be sure the RNA quantitation using UV absorbance is performed with 10 mM Tris-HCl, pH 7.0 (page 21) to accurately measure the UV absorbance.	
	Loss of pellet during to ethanol precipitation	 Remove supernatant from RNA pellet carefully. Use RiboMinus[™] Concentration Module to concentrate the 	
		RiboMinus™ RNA.	
Incomplete removal of rRNA	Too much total RNA used	The protocols in this manual are designed to purify RiboMinus™ RNA from 2-10 µg total RNA. If you are using more than 10 µg total RNA, divide the sample into two sample aliquots, each containing <10 µg total RNA for RiboMinus™ RNA purification.	
	Low amount of magnetic beads or probe used	Be sure to use the recommended amounts of RiboMinus™ Probe and RiboMinus™ Magnetic Beads for efficient removal of rRNA.	

Troubleshooting, Continued

Problem	Cause	Solution	
Incomplete removal of rRNA	Improper handling or drying of beads	To obtain the best results with RiboMinus™ Magnetic Beads, follow the guidelines on page 12 for washing and mixing the beads, and aspirating the supernatant. Do not allow the beads to dry as drying reduces the bead efficiency.	
RNA degraded	RNA contaminated with RNase	Follow the guidelines on page 7 to prevent RNase contamination.	
	Poor quality starting materials	Always use fresh samples or samples frozen at -80°C for isolation of total RNA. Be sure to check the quality of your total RNA prior to use.	
Genomic DNA contamination	Total RNA contained genomic DNA	Perform DNase I digestion with the total RNA sample to remove any genomic DNA contamination before performing RiboMinus™ RNA purification.	
Inhibition of downstream enzymatic reactions	Presence of ethanol in purified RNA sample	Traces of ethanol from the RiboMinus™ Concentration Module Wash Buffer (W5) or from ethanol precipitation can inhibit downstream enzymatic reactions.	
		To remove Wash Buffer (W5), discard Wash Buffer flow through from the collection tube. Reinsert the spin column into the collection tube and centrifuge the spin column at maximum speed for 2 minutes to completely dry the column.	
		 If performing ethanol precipitation, make sure that ethanol is evaporated before resuspending the RiboMinus™ RNA pellet in RNase- free water. 	

Technical Service

World Wide Web



Visit the Invitrogen Web Resource using your World Wide Web browser. At the site, you can:

- Get the scoop on our hot new products and special product offers
- View and download vector maps and sequences
- Download manuals in Adobe® Acrobat® (PDF) format
- Explore our catalog with full color graphics
- Obtain citations for Invitrogen products
- Request catalog and product literature

Once connected to the Internet, launch your Web browser (Internet Explorer 5.0 or newer or Netscape 4.0 or newer), then enter the following location (or URL):

http://www.invitrogen.com

...and the program will connect directly. Click on underlined text or outlined graphics to explore. Don't forget to put a bookmark at our site for easy reference!

MSDS Requests

To request an MSDS, visit our Web site at www.invitrogen.com. On the home page, go to 'Technical Resources', select 'MSDS', and follow instructions on the page.

Purchaser Notification

Limited Warranty

Invitrogen is committed to providing our customers with highquality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, please contact our Technical Service Representatives. Invitrogen warrants that all of its products will perform according to the specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's <u>liability</u> only to the cost of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives. Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.

References

- Braasch, D. A., and Corey, D. R. (2001). Locked Nucleic Acid (LNA): Fine-tuning the Recognition of DNA and RNA. Chem Biol. 1, 1-7.
- McTigue, P. M., Peterson, R. J., and Kahn, J. D. (2004). Sequence-dependent Thermodynamic Parameters for Locked Nucleic Acid (LNA)-DNA Duplex Formation. Biochemistry. 43, 5388-5405.
- Ruan, Y., Le Ber, P., Ng, H., and Liu, E. (2004). Interrogating the Transcriptome. Trends Biotechnol. 22, 23-30.

©2005 Invitrogen Corporation. All rights reserved.

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

RNase $AWAY^{\otimes}$ and TRIzol $^{\otimes}$ are registered trademarks of Molecular Bio-Products, Inc. LNA $^{\bowtie}$ is a trademark of Exiqon A/S. LabChip $^{\otimes}$ is a registered trademark of Caliper Life Sciences, Inc.



Corporate Headquarters:

Invitrogen Corporation 1600 Faraday Avenue Carlsbad, California 92008 Tel: 1 760 603 7200 Tel (Toll Free): 1 800 955 6288 Fax: 1 760 603 7229

Email: tech_service@invitrogen.com

European Headquarters:

Invitrogen Ltd 3 Fountain Drive Inchinnan Business Park Paisley PA4 9RF, UK Tel (Free Phone Orders): 0800 269 210 Tel (General Enquiries): 0800 5345 5345 Fax: +44 (0) 141 814 6287 Email: eurotech@invitrogen.com

International Offices:

Argentina 5411 4556 0844 Australia 1 800 331 627 Austria 0800 20 1087 Belgium 0800 14894 Brazil 0800 11 0575 Canada 800 263 6236 China 10 6849 2578 Denmark 80 30 17 40

France 0800 23 20 79
Germany 0800 083 0902
Hong Kong 2407 8450
India 11 577 3282
Italy 02 98 22 201
Japan 03 3663 7974
The Netherlands 0800 099 3310
New Zealand 0800 600 200
Norway 00800 5456 5456

Spain & Portugal 900 181 461 Sweden 020 26 34 52 Switzerland 0800 848 800 Taiwan 2 2651 6156 UK 0800 838 380 For other countries see our Web site

www.invitrogen.com

