USER GUIDE



Dynabeads[®] mRNA DIRECT[™] Kit

For the isolation of pure mRNA directly from crude samples

Catalog numbers 61011 and 61012

Revision Date June 2012 (Rev. 009)



For Research Use Only. Not for human or animal therapeutic or diagnostic use.

Contents

Kit Contents and Storage	v
Description of the System	1
About the Kit	1
Experimental Overview	3
Methods	4
General Guidelines	4
Sample Preparation	5
Prepare Dynabeads [®] Oligo (dT) ₂₅	8
Direct mRNA Isolation Protocol	9
Large Scale Isolations	11
Elimination of rRNA Contamination	12
Troubleshooting	
Appendix	16
Accessory Products	16
Sample Types from which mRNA has been Isolated Using Dynabeads [®] Oligo (dT) ₂₅	17
Technical Support	
Purchaser Notification	
References	

Kit Contents and Storage

Storage

Upon receipt, store all components of the Dynabeads[®] mRNA DIRECT[™] Kit at 2°C to 8°C. **Do not freeze** the Dynabeads[®] Oligo (dT)₂₅.

Keep Dynabeads[®] in liquid suspension during storage and all handling steps, as drying will result in reduced performance. Resuspend thoroughly before use.

Precautions should be taken to ensure that RNase or microbial contamination of the kit components does not occur.

This product contains 0.02% sodium azide as a preservative.



Warning! GENERAL CHEMICAL HANDLING.

For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Sodium azide may react with lead and copper plumbing to form highly explosive metal azides.

Kit Contents and Storage, continued

Kit contents The components included in the Dynabeads[®] mRNA DIRECT[™] Kit are listed in the following table. The suspension of Dynabeads[®] Oligo (dT)₂₅ and the buffers provided are produced and packed under RNase-free conditions. All kit reagents are of analytical grade and are RNase-free.

Sufficient reagents are provided in the kit to perform 20 (Cat. no. 61011) or 40 (Cat. no. 61012) isolations.

Component	Cat. no. 61011	Cat. no. 61012
Dynabeads[®] Oligo (dT) 25 [*] (≈5 mg/mL, supplied in PBS pH 7.4)	5 mL	10 mL
Lysis/Binding Buffer 100 mM Tris-HCl, pH 7.5 500 mM LiCl 10 mM EDTA, pH 8 1% LiDS 5 mM dithiothreitol (DTT)	30 mL	60 mL
Washing Buffer A 10 mM Tris-HCl, pH 7.5 0.15 M LiCl 1 mM EDTA 0.1% LiDS	60 mL	120 mL
Washing Buffer B 10 mM Tris-HCl, pH 7.5 0.15 M LiCl 1 mM EDTA	30 mL	60 mL
10 mM Tris-HCl pH 7.5 (Elution Buffer) *Contains 0.02% sodium azide as preservative.	15 mL	15 mL

*Contains 0.02% sodium azide as preservative.

Product Use

For Research Use Only. Not for human or animal therapeutic or diagnostic use.

Description of the System

About the Kit

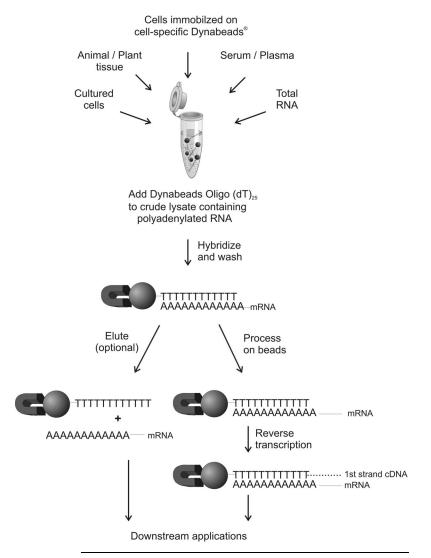
System The isolation protocol relies on base pairing between	Introduction	This product is designed for simple and rapid isolation of pure, intact polyadenylated (polyA) mRNA directly from the crude lysate of animal and plant cells and tissues. The isolated mRNA is suitable for use in all downstream applications.
overviewthe polyA residues at the 3' end of most mRNA, and the oligo (dT)25 residues covalently coupled to the surface of the Dynabeads®. Other RNA species lacking a polyA tail will not hybridize to the beads and are readily washed away. RNase inhibiting agents in the Lysis/Binding Buffer together with stringent hybridization and washing conditions ensure the isolation of pure, intact mRNA from crude samples rich in RNase, without the use of strong chaotropic agents. The protocol is flexible and can easily be scaled up or down to suit all sample sizes. It has successfully been used in the isolation of mRNA from single cells (Karrer, et al., 1995). The high capture efficiency facilitates detection of mRNA by reverse transcriptase (RT)-PCR from highly specialized cells (e.g., isolated 	System overview	the polyA residues at the 3' end of most mRNA, and the oligo (dT) ₂₅ residues covalently coupled to the surface of the Dynabeads [®] . Other RNA species lacking a polyA tail will not hybridize to the beads and are readily washed away. RNase inhibiting agents in the Lysis/Binding Buffer together with stringent hybridization and washing conditions ensure the isolation of pure, intact mRNA from crude samples rich in RNase, without the use of strong chaotropic agents. The protocol is flexible and can easily be scaled up or down to suit all sample sizes. It has successfully been used in the isolation of mRNA from single cells (Karrer, et al., 1995). The high capture efficiency facilitates detection of mRNA by reverse transcriptase (RT)-PCR from highly specialized cells (e.g., isolated from a heterogeneous sample by immunomagnetic separation). In addition, the protocol has been successfully used to isolate mRNA from a wide variety of tissues of mammalian, fish, amphibian,

About the Kit, continued

System overview, continued	For many applications elution of the mRNA from the beads is not required as the beads do not interfere with downstream enzymatic reactions. The beadbound oligo (dT) ₂₅ can also function as a primer for RT and synthesis of first-strand cDNA (ref. 60), allowing the construction of solid-phase cDNA libraries and solid-phase RT-PCR.		
Binding Capacity	1 mg of Dynabeads [®] Oligo $(dT)_{25}$ beads (200 µL) binds up to 2 µg of mRNA. A typical mammalian cell contains about 10–30 pg of total RNA, from which 1–5% is mRNA.		
Expected Yield	The information in the following table is intended as a rough guide to the expected total RNA content of selected tissues, as well as appropriate bead and buffer volumes.		
	Cell Types & Quantity Estimated Total RNA content (1–5% is mRNA)		
	Cell Types & Quantity		
	Cell Types & Quantity Single Mammalian Cell	Total RNA content (1–5% is mRNA)	
		Total RNA content	
	Single Mammalian Cell	Total RNA content (1–5% is mRNA) 10–30 pg	
	Single Mammalian Cell 50 mg of muscle tissue	Total RNA content (1–5% is mRNA) 10–30 pg 50–80 μg	
	Single Mammalian Cell 50 mg of muscle tissue 50 mg of liver tissue	Total RNA content (1–5% is mRNA) 10–30 pg 50–80 μg 400 μg	

Experimental Overview

mRNA Isolation Workflow Outline of the protocol for isolating mRNA from a crude starting sample using Dynabeads[®] Oligo(dT)₂₅. The isolated mRNA is suitable for use in all downstream molecular biology applications.



Methods

General Guidelines

General Guidelines

- Keep Dynabeads[®] Oligo (dT)₂₅ in liquid suspension during storage and all handling steps. Resuspend thoroughly before use.
- Work RNase free and wear gloves.
- Precipitate may form in the buffers. Dissolve precipitate before use by warming to room temperature and mixing thoroughly.
- Bring Lysis/Binding Buffer and Washing Buffers A and B to room temperature prior to use.
- 10mM Tris-HCl pH 7.5 (Elution Buffer) should be stored on ice or at 2°C to 8°C prior to use.
- Thorough resuspension of the beads/mRNA complex during washing and complete removal of the washing buffer at each step will prevent carryover of LiDS and other salts to the downstream reaction. Transferring the beads/ mRNA complex to new tubes before the last washing step will further reduce LiDS carry over. LiDS is a strong inhibitor of enzymatic reactions.

Sample Preparation

Introduction	mRNA content of cells and tissues vary depending on the source of the material and RNA expression levels at the time of tissue/cell harvest. Dynabeads [®] mRNA DIRECT [™] Kit protocols can be scaled up or down to suit specific sample source and quantity.
Required	Components required but not supplied:
Materials	 RNase-free micropipettors, disposable pipette tips (aerosol resistant preferred) and micro- centrifuge tubes
	 DynaMag[™]-2 magnet, see Accessory Products (page 16) for ordering information or visit www.lifetechnologies.com/magnets for magnet recommendations
	• Mixer allowing both tilting and rotation
	• Water bath or heating block (for optional elution)
	• Microcentrifuge capable of achieving $>10,000 \times g$
	Components required for isolation of mRNA from tissue samples:
	Liquid Nitrogen
	Mechanical or manual tissue grinder
	• Syringe and 21 gauge needle
	Components supplied with the kit:
	 Dynabeads[®] Oligo (dT)₂₅
	Lysis/Binding Buffer
	Washing Buffer A
	Washing Buffer B
	• Elution Buffer (10 mM Tris-HCl, pH 7.5)
	Continued on next page

Sample Preparation, continued

Solid Plant or Animal Tissue

1. Aliquot (weigh) the animal or plant tissue while frozen, to avoid mRNA degradation. Ideally the tissue should be weighed and aliquoted before freezing. Do not exceed the specified amount of tissue (see the following table), as using too much tissue reduces the mRNA yield and purity.

Component	Std (mg)	Micro (mg)	Mini (mg)	Maxi (mg)
Plant Tissue	20–100	≤4	4–20	100-400
Animal Tissue	10–50	≤2	2–10	50-200

- 2. Grind frozen tissue in liquid nitrogen. Work quickly.
- Transfer the frozen powder to the appropriate volume of Lysis/Binding Buffer and homogenize until complete lysis is obtained (≈1–2 min). Rapid lysis in the Lysis/Binding Buffer is critical for preventing mRNA degradation.

Std	Micro	Mini	Maxi
1250 µL	300 µL	300 µL	5 mL

4. A DNA-shear step is recommended for samples containing over 5×10^5 cells.

Force the lysate through a 21 gauge needle 3–5 times using a 1–2 mL syringe to shear the DNA. The reduction in viscosity should be noticeable.

Note: Repeated shearing causes foaming of the lysate due to detergent in the buffer, however, this should not affect the mRNA yield. The foam can be reduced by a 30 second centrifugation. The lysate can be frozen and stored at –80°C for later use.

Sample Preparation, continued

Cultured Cells or Cell Suspension

1. Pellet the appropriate number of cells (indicated in the following table), by centrifugation (e.g., $400 \times g$ for 8 minutes at 4°C).

Std	Micro	Mini	Maxi
$1-4 \times 10^6$	≤150,000	$0.15 - 1 \times 10^{6}$	$4-20 \times 10^{6}$

- 2. Wash the pellet by resuspending in phosphatebuffered saline (PBS). Pellet cells by centrifugation again. Use the cell pellet immediately, or freeze in liquid nitrogen or at -80°C for later use.
- 3. Add the appropriate volume of Lysis/Binding Buffer to either a frozen cell pellet or to a fresh cell pellet.

Std	Micro	Mini	Maxi
1250 µL	300 µL	300 µL	5 mL

- 4. Perform a repeated passage of the solution through a pipette tip to obtain complete lysis. The release of DNA during lysis results in a viscous solution which confirms complete lysis.
- 5. A DNA-shear step is recommended for samples containing over 5×10^5 cells.

Force the lysate through a 21 gauge needle 3–5 times using a 1–2 mL syringe to shear the DNA. The reduction in viscosity should be noticeable.

Note: Repeated shearing causes foaming of the lysate due to detergent in the buffer, however, this should not affect the mRNA yield. The foam can be reduced by a 30 second centrifugation. The lysate can be frozen and stored at –80°C for later use.

Prepare Dynabeads® Oligo (dT)25

Prepare Dynabeads® Oligo (dT)25

- 1. Resuspend Dynabeads[®] Oligo(dT)₂₅ thoroughly before use.
- Transfer the desired volume of beads from the stock tube to an RNase-free 1.5-mL microcentrifuge tube and place the tube on a magnet (e.g., DynaMag[™]-2).

Std	Micro	Mini	Maxi
250 µL	10 µL	50 µL	1 mL

- 3. After 30 seconds (or when the suspension is clear), remove the supernatant.
- 4. Remove the tube from the magnet and wash the beads by resuspending in an equivalent volume of fresh Lysis/Binding Buffer.

Optional: For very small bead volumes (mini- and micro- isolations) use 50–100 µL wash volume to ease handling.

5. Proceed to Direct mRNA Isolation Protocol, page 9.

Direct mRNA Isolation Protocol

Direct mRNA Isolation Protocol

- Remove the Lysis/Binding Buffer from the prewashed Dynabeads[®] Oligo(dT)₂₅ (Prepare Dynabeads[®] Oligo (dT)₂₅, step 4) by placing on the magnet for 30 seconds, or until the suspension is clear and discard the supernatant.
- 2. Remove the microcentrifuge tube from the magnet and add the appropriate volume of sample lysate indicted in the following table.

Std	Micro	Mini	Maxi
1250 µL	300 µL	300 µL	5 mL

- 3. Pipet to resuspend the beads completely in the sample lysate.
- Incubate with continuous mixing (rotating or roller mixer) for 3–5 min at room temperature to allow the polyA tail of the mRNA to hybridize to the oligo (dT)₂₅ on the beads. Note: Increase the incubation time if the solution is viscous.
- 5. Place the vial on the magnet for 2 minutes and remove the supernatant. If the solution is noticeably viscous, increase the time to approximately 10 minutes.
- 6. Wash the beads/mRNA complex two times with the appropriate volume of Washing Buffer A (see the following table) at room temperature. Use the magnet to separate the beads from the solution between each washing step.

Std	l	Micro	Mini	Maxi
1–2 mI	,	600 µL	600 µL	10 mL

Direct mRNA Isolation Protocol, continued

Direct mRNA Isolation Protocol, continued

7.

Wash the beads/mRNA complex once with the appropriate volume of Washing Buffer B (see the following table) at room temperature. Use the magnet to separate the beads from the solution.

Std	Micro	Mini	Maxi
1–1.5 mL	300 µL	300 µL	5 mL

8. If the isolated mRNA is to be used in enzymatic downstream applications (e.g., RT-PCR), one extra wash in Washing Buffer B is recommended. This should be followed by a final wash in the enzymatic buffer to be used (e.g., RT-PCR buffer without the enzyme or primers).

Note: Perform cDNA synthesis as recommended by the manufacturer of the reverse transcriptase. When using a thermostable reverse transcriptase and the bead-bound oligo (dT) as primer for first strand cDNA synthesis, an initial incubation at 50°C for 5 minutes is necessary before proceeding at the recommended temperature.

9. If elution of mRNA from the beads is desired, add an appropriate volume (see the following table) of 10 mM Tris-HCl pH 7.5 (Elution Buffer) and incubate at 65°C to 80°C for 2 minutes. Immediately place the tube on the magnet, transfer the supernatant containing the mRNA to a new RNase-free tube and place this tube on ice.

Std	Micro	Mini	Maxi
10–25 μL	10 µL	10 µL	50–100 μL

Large Scale Isolations

Reuse of Dynabeads® Oligo (dT)25	Note: The buffers supplied with the kit may not be sufficient for large scale mRNA isolations. Multiple isolations from the same sample can be performed by reusing Dynabeads [®] Oligo(dT) ₂₅ after mRNA elution. Simply follow the protocol described in Direct mRNA Isolation Protocol (page 9).
	 After elution of the mRNA, wash the beads once in Lysis/Binding Buffer (Prepare Dynabeads[®] Oligo (dT)₂₅, page 8).
	• Add a new lysate sample to the beads and continue the isolation as usual. Alternatively, washed beads can be re-applied to the same sample lysate until all the mRNA has been captured.

Elimination of rRNA Contamination

Elimination of rRNA Contamination	bee apj tra int	some cases trace amounts of ribosomal RNA have en observed in the mRNA samples. For many plications such as Northern blotting and RT-PCR, ce amounts of rRNA contamination will not erfere with the analysis or interpretation of the sults.	
	However, for other applications such as cDNA library construction and microarray analysis, rRNA contamination should be avoided.		
	ext sar iso wa	posomal RNA is effectively eliminated by re- cracting the mRNA from the eluate. Reuse of the me Dynabeads [®] Oligo (dT) ₂₅ used for the original lation is recommended. If new beads are used, ush the beads in 50 mM sodium pyrophosphate fore the isolation of mRNA.	
	1.	Follow the Direct mRNA Isolation Protocol (page 9). Elute the mRNA in Elution Buffer.	
	2.	Transfer the eluted mRNA to a new RNase-free tube and place on ice. Do not discard the beads.	
	3.	Wash the beads two times in Washing Buffer B (Direct mRNA Isolation Protocol , step 6 on page 10).	
	4.	Dilute the eluted mRNA with 4 times its volume of Lysis/Binding Buffer (e.g., if the mRNA is eluted in 20 μ L, add 80 μ L of Lysis/Binding Buffer).	
	5.	Remove the Washing Buffer B from the beads, by placing the tube on the magnet, and add the diluted mRNA.	
	6.	Incubate with mixing at room temperature for 3–5 minutes.	
	7.	Continue with the Direct mRNA Isolation Protocol (page 9) starting at step 4.	
		(I0)	

Observation	Cause	Solution
Clumping of beads during	DNA in the sample lysate has not been	Pipet the solution several times through a 1-mL pipette.
incubation with sample lysate.	completely sheared.	Increase force/number of passages through the needle in future shearing steps.
mRNA is contaminated with DNA.	Incomplete DNA shearing.	Increase the force and/or the number of passages through the needle in the DNA shearing step.
	Incomplete removal of sample lysate after hybridization step, and subsequent carry over to wash and elution steps.	Completely remove the sample lysate after hybridization.
	Inefficient washing.	Make sure the beads/mRNA complex is fully resuspended in washing buffer.
	Incomplete removal of wash buffers.	Completely remove the sample/washing buffers.
	Sample-to-beads ratio too high.	Dilute sample lysate or increase the amount of beads.
		Re-extract the mRNA from the eluate.
mRNA yield lower than expected.	Inefficient elution of mRNA from the beads.	Increase the elution volume/time/ temperature or perform the elution step two times, pooling the eluate.
	Beads-to-sample ratio is too low.	Increase the amount of beads.
	Cells/tissue incompletely lysed.	Repeat the homogenization step.

Troubleshooting, continued

Observation	Cause	Solution
The beads- cDNA complex is clumped and sticking to the tubes after reverse transcription.	Non-specific electrostatic interactions between the cDNA molecules and the plastic materials of the tubes/pipette tips.	 Add BSA (0.2–1.0% final concentration) to the reverse transcription mix before performing the cDNA synthesis. This is to reduce clumping of the beads for a more efficient cDNA synthesis. Note: Use best possible BSA quality. Where appropriate, add 0.05% Tween-20 to the reaction buffers. Alternatively, Dilute the beads/cDNA solution (after reverse transcription) with an equal volume of the 1X reverse transcription reaction buffer containing 0.05% Tween-20. Mix by pipetting and transfer the suspension to a new tube. If there are any remaining beads stuck to the tube walls, remove by washing with a fresh aliquot of buffer containing Tween-20. Pool these beads with the original bead suspension. Place the pooled beads on a magnet and remove the supernatant. Wash 2–3 times with the buffer containing Tween-20. Store the solid-phase cDNA library in an appropriate buffer containing 0.05% Tween-20.

Troubleshooting, continued

Observation	Cause	Solution
Unable to detect specific	The beads-to- sample ratio is too low.	Increase the amount of beads.
mRNA molecules.	Inappropriate sample volume.	Reduce sample volume/increase sample concentration.
	Hybridization time too short.	Increase hybridization incubation time to 10–15 min.

Appendix

Accessory Products

IntroductionThe following products may be used with the
Dynabeads[®] mRNA DIRECT[™] Kit. For details, visit
www.lifetechnologies.com or contact Technical
Support (see page 21).

Product	Quantity	Cat. no.
Dynabeads [®] Oligo(dT) ₂₅	5 mL	61005
DynaMag [™] -2 Magnet	Each	12321D
SampleRack (for DynaMag [™] -2)	Each	12322D
DynaMag [™] -15 Magnet	Each	12301D
DynaMag [™] -96 Side Skirted	Each	12027
HulaMixer [®] Sample Mixer	Each	15920D
SuperScript [®] III Reverse Transcriptase	2000 units	18080-093
SuperScript [®] III First-Strand Synthesis System	50 reactions	18080-051
SuperScript [®] III One-Step RT-PCR System with Platinum [®] Taq	100 reactions	12574-035
SuperScript [®] VILO [™] cDNA Synthesis Kit	50 reactions	11754-050
UltraPure [™] DNase/RNase-free Distilled Water	500 mL	10977-015

Sample Types from which mRNA has been Isolated Using Dynabeads® Oligo (dT)₂₅

Tissue	Species	References
Adrenals	Rat	4
Brain	Mouse, Trout	12, 4
Brain (cerebral cortex, preoptic area, dentate gyrus)	Rat	13, 14, 15
Cartilage	Human	16
Organ of Corti and spiral ganglion	Guinea Pig, Rat	17
Ear (cochleae)	Mouse	11
Eggs	Trout	4
Gut (paraffin embedded)	Human	18
Heart	Rat	13, 19*
Hypothalamus	Rat	15
Kidney	Rat	13
Kidney (glomerular preparations)	Human	20
Liver (paraffin embedded)	Human	18
Liver	Rat, trout, Xenopus	13, 19*, 4
Lung (paraffin embedded)	Human	18
Lung	Rat	13
Muscle	Rat, trout	19*, 4
Nematode (frozen rehydrated cysts)	Globodera rostochiensis	58
Ovaries	Trout, Xenopus	4
Pancreas	Rat	19*
Paraffin embedded lung, liver, gut	Human	18
Paraffin embedded keratinocytes	Human	21
Pituitary	Rat	22
Plasma	Human	59
Pronephos	Trout	4
Skin (dried)	Frog	57
Spleen	Rat	4, 13
Trematode	Schistosoma mansoni	23
Whole insect	Drosophila	4

Animal Tissues Isolated mRNA from animal tissues

*Lysis buffer with 4 M urea and 1 % SDS

Sample Types from which mRNA has been Isolated Using Dynabeads® Oligo (dT)25, continued

Tissue	Species	References
Whole plants	Arabidopsis thaliana, Rice, Oryza sativa	23, 24, 25, 26, 27
Bud	Tobacco	28
Epidermal leaf cell (single cells)	Tomato	1
Embryos	Maize	29
Flowers	Maize, Tobacco	30
Guard cell in leaf (single cells)	Tomato	1
Leaves	Barley, <i>Brassica oleracea</i> , Maize, Potato, Tobacco, Tomato	31, 4, 32, 33, 29, 27, 34, 28, 1
Ovules	Maize	29
Roots	Barley, <i>Brassica oleracea,</i> Spruce, Maize	31, 33, 6, 29
Seed aleurone	Barley	31, 4, 35, 36, 32
Seed endosperm	Barley	35, 36, 32
Seed embryos	Barley	31, 4, 35, 36, 32
Seedlings	Maize, Tobacco	29, 28
Single leaf cells	Tomato	1
Stem	Tobacco	28
Stigma	Brassica oleracea	33, 37
Stolon tips	Potato	34

Plant Tissues Isolated mRNA from plant tissues

Sample Types from which mRNA has been Isolated Using Dynabeads® Oligo (dT)25, continued

Cells

Isolated mRNA from different types of cells

Cell type/cell line	Species	References
Chondrocytes	Human	16
Cervical cancer cells (HeLa)	Human	38, 39
Colon carcinoma cell line (COLO320)	Human	40
Fibroblast cells line (ST-1 and SKB-1)	Human	41, 9
Fibroblast (D551)	Human	8,9
Fibroblast (RTG-2)	Trout	4
Endothelial cells (umbilical cord)	Human	8,9
Hepatocyte cell line (HepG2)	Human	8, 39
Keratinocytes	Human	42, 21
Langerhans cells	Human	42
Lymphoblast B-cell lines (Reh, Daudi, HL-60, IM9)	Human	8,4,39,43
Mamma carcinoma cells (MCF7)	Human	38, 39
Mamma carcinoma (T47D)	Human	40
Monocytes	Human	44
Pancreas, insulinoma Rinm5F cells	Rat	45
Peripheral blood mononuclear cells (PBMC)	Human	46
Peritoneal exuduate cells	Human	42
Placental cell line (AMA)	Human	38, 39
T-cells/T-cell clones	Human	2, 47, 48, 49, 50
Yeast (Saccharomyces cerevisiae, Hansenula polymorpha)	In soil samples	51
Yeast (Saccharomyces cerevisiae)	Culture	52

Sample Types from which mRNA has been Isolated Using Dynabeads® Oligo (dT)25, continued

Viruses

Direct isolation of viral polyA RNA with Dynabeads $^{\ensuremath{\$}}$ Oligo (dT)₂₅

Starting Material	Virus	References
Cells in bronchoalveolar washes	HIV-1	47*
Cerebrospinal fluid	HIV-1	53*
Cell line	HTLV-I/II	54**
Peripheral blood mononuclear cells (PBMC)	HIV-1	47*
Plasma	HIV-1/HIV-2	54*,**
Serum	HIV-1	55*, 56*, 54*, **
T-lymphocytes cell line (CD4+)	HIV-1	47*

*Lysis/binding buffer: 1 M LiCl, 2% SDS, 2XTE, 50 µg tRNA, Vanadyl ribonucleosyl complexes.

**Lysis/binding buffer: 4 M GTC, 0.5% sarkosyl, 1% DTT, 0.5 M LiCl, 0.1 M Tris pH8.

Technical Support

Obtaining Support	 For the latest services and support information for all locations, go to www.lifetechnologies.com. At the website, you can: Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities Search through frequently asked questions (FAQs) Submit a question directly to Technical Support (techsupport@lifetechnologies.com) Search for user documents, Safety Data Sheets (SDSs), vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents Obtain information about customer training Download software updates and patches
Safety Data Sheets (SDS)	Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/support .
Certificate of Analysis	The Certificate of Analysis is available at www.lifetechnologies.com/support . Search for the Certificate of Analysis by product lot number, which is printed on the box.
Limited Product Warranty	Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions . If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support .

Purchaser Notification

Limited Use Label License: Research Use Only	The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.
General	Life Technologies AS complies with the Quality
Information	System Standards ISO 9001:2008 and ISO 13485:2003.

References

- Karrer EE, et al. (1995) In situ isolation of mRNA from individual plant cells: Creation of cell-specific cDNA libraries. Proc Natl Acad Sci USA 92:3814-3818.
- Raineri I, et al. (1991) Improved efficiency for single-sided PCR by creating a reusable pool of first strand cDNA coupled to a solid phase. Nucleic Acids Research 19(14):4010.
- Lambert KN and Williamson VM. (1993) cDNA library construction from small amounts of RNA using paramagnetic beads and PCR. Nucleic Acids Research 21(3):775-776.
- Jakobsen KS, et al. (1994) Direct mRNA isolation using Magnetic Oligo(dT) Beads: A protocol for all types of cell cultures, animal and plant tissues. Advances in Biomagnetic Separation, Ed. Uhlèn M, Hornes E and Olsvik Ø, Eaton Publishing 61-71.
- 5. Rodriguez IR and Chader GJ. (1992) A novel method for the isolation of tissuespecific genes. Nucleic Acids Research 20(13):3528.
- 6. Sharma P, et al. (1993) PCR-based construction of subtractive cDNA library using magnetic beads. BioTechniques 15(4):610-611.
- 7. Schraml P, et al. (1993) cDNA subtraction library construction using a magnetassisted subtraction technique (MAST). Trends in Genetics 9(3):70-71.
- Aasheim HC, et al. (1994) A simple subtraction method for the isolation of cellspecific genes using magnetic monodisperse polymer particles. Biotechniques 16(4):716-721.
- 9. Coche T, et al. (1994) Generation of an unlimited supply of a subtracted probe using magnetic beads and PCR. Nucleic Acids Research 22(7):1322-1323.
- 10. Borgnes A, et al. Detection of isolated colon carcinoma cells in peripheral blood- or bone marrow mononuclear cell suspensions. Micrometastasis Meeting Munich, June 1996 (Abstract).
- 11. Dazert SB, et al. (2007) Hearing development and spiral ganglion neurite growth in VASP deficient mice. Brain Research 1178:73-82.
- 12. Moyal M, et al. (1992) Mutations in the UL53 gene of SV-1 abolish virus neurovirulence to mice by the intracerebral route of infection. Virus Research 26:99-112.
- 13. Pines G, et al. (1992) Cloning and expression of a rat brain L- glutamate transporter. Nature 360:464-467.
- 14. Alme MN, et al. (2007) Chronic Fluoxetine Treatment Induces Brain Region-Specific Upregulation of Genes Associated with BDNF-Induced Long-Term Potentiation. Neuronal Plasticity:1-9.
- 15. Kim K, et al. (1994) Competitive PCR for quantification of gonadotropinreleasing hormone mRNA level in a single micropunch of the rat preoptic area. Mol Cel Endocrinol 97:153-158.

- 16. Winterpacht A, et al. (1994) Alter-native splicing as the result of a type II procollagen gene (COL2A1) mutation in a patient with Kneist Dysplasia. Hum Mol Genet 3(10):1891-1893.
- 17. Safieddine S and Wenthold RJ. (1997) The glutamate receptor subunit d1 is highly expressed in hair cells of the auditory and vestibular systems. J Neurosci 17:7523-7531.
- 18. de Andrès B, et al. (1995) Improved method for mRNA extraction from paraffin-embedded tissues. BioTechniques 18(1):42-44.
- 19. Hengerer B. (1993) A rapid procedure for mRNA extraction from a large number of samples. Biotechniques 14(4):522-524.
- Bicknell G, et al. (1996) Separate extraction of mRNA from glomerular epithelial and mesangial cells. Pathological Society of Great Britain and Ireland, 172nd Meeting.
- 21. Shamsher MK, et al. (1995) Novel mutations in keratin 16 gene underly focal non-epidermolytic palmoplantar keratoderma (NEPPK) in two families. Hum Mol Genet 4(10):1875-1881.
- Schauder B, et al. (1994) Cloning of a cDNA encoding an ectoenzyme that degrades thyrotropin-releasing hormone. Proc Natl Acad Sci USA 91:9534-9538.
- 23. Schüssler P, et al. (1995) Combined isolation of nucleic acids and protein from small amounts of tissue. Trends in Genetics 11(10):378-379.
- 24. Rubin D. Magnetic Bead Isolation of mRNA from Plant Tissue. In: The Red Book Bulletin, Current Protocols in Molecular Biology, Subscriber's Notebook, Anonymous 1992, p. 1, Suppl. 20.
- 25. Ivanov R. et al. (2008) EFFECTOR OF TRANSCRIPTION2 is involved in xylem differentiation and includes a functional DNA single strand cutting domain. Developmental Biology (313):93-106.
- 26. Magrelli A, et al. (1994) Splicing of the rolA transcript of Agrobacterium rhizogenes in Arabidopsis. Science 266:1986-1988.
- Zhu QH, et al. (2007) compact shoot and leafy head 1, a mutation affects leaf initiation and developmental transition in rice (Oryza sativa L.). Plant Cell Reports (26):421-427.
- 28. Hewelt A, et al. (1994) Promoter tagging with a promoterless ipt gene leads to cytokinin-induced phenotypic variability in transgenic tobacco plants: implications of gene dosage effects. Plant Journal 6(6):879- 891.
- 29. Breton C, et al. (1995) PCR-generated cDNA library of transition-stage maize embryos: cloning and expression of calmodulin genes during early embryogenesis. Plant Mol.Biol. 27:105-113.

- 30. Allen RL and Lonsdale DM. (1993) Molecular characterization of one of the maize poly-galacturonase gene family members which are expressed during late pollen development. Plant Journal 3(2):261- 271.
- Jakobsen KS, et al. (1990) Purification of mRNA directly from crude plant tissues in 15 minutes using oligo dT microspheres. Nucleic Acids Research 18(12):3669.
- 32. Stacy RAP, et al. (1995) Evolution of the group 1 late embryogenesis abundant (Lea) genes: analysis of the Lea B19 gene family in Barley. Plant Mol Biol 28:1039-1054.
- Kumar V and Trick M. (1994) Expression of the S-locus receptor kinase multigene family in Brassica oleracea. Plant Journal 6(6):807-813.
- Bachem CWB, et al. (1994) Antisense expression of polyphenol oxidase genes inhibits enzymatic browning in potato tubers. Biotechnology 12:1101-1105.
- 35. Espelund M, et al. (1992) Late embryogenesis-abundant genes encoding proteins with different numbers of hydrophilic repeats are regulated differentially by abscisic acid and osmotic stress. Plant Journal 2(2):241-252.
- 36. Aalen R, et al. (1994) Transcripts encoding oleosin and a dormancyrelated protein are present in both the aleurone layer and the embryo of developing barley (Hordeum vulgare L.) seeds. Plant Journal 5(3):385-396.
- 37. Giranton JL, et al. (1995) The S locus receptor kinase gene encodes a soluble glycoprotein corresponding to the SRK extracellular domain in Brassica oleracea. Plant Journal 8(6):827-834.
- Larsen F, et al. (1993) A tight cluster of five unrelated human genes on chromosome 16q22.1. Hum Mol Genet 2(10):1589-1595.
- 39. Luna L, et al. (1994) Molecular cloning of a putative novel human bZIP transcription factor on chromosome 17q22. Genomics 22:553-562.
- 40. Maas RA, et al. (1995) Immuno-magnetic purification of human breast carcinoma cells allows tumor-specific detection of multidrug resistance gene-1 mRNA by reverse transcriptase polymerase chain reaction in fine needle aspirates. Lab Invest 72(6):760-764.
- 41. Slupphaug G, et al. (1991) Cell cycle regulation and in vitro hybrid arrest analysis of the major human uracil-DNA glycosylase. Nucleic Acids Research 19:5131-5137.
- 42. Sprecher E and Becker Y. (1992) Detection of IL-1, TNF- and IL-6 gene transcription by the polymerase chain reaction in keratinocytes: Langerhans cells and peritoneal exudate cells during infection with herpes simplex virus-1. Arch Virology 126:253-269.

- 43. Day S and Ross J. (1995) mRNA based method for typing the HLA-C locus gene by automated sequencing based typing. Eur J Immunogenet 23(1):92.
- 44. Osnes LTN, et al. (1995) Lipopolysaccharide activation of human monocytes mediated by CD14, results in a coordinated synthesis of tissue factor TNFalpha and IL-6. J Endotoxin Res 2:27-35.
- 45. Bristulf J, et al. (1994) Interleukin-1 stimulates the expression of type I and type II interleukin-1 receptors in the rat insulinoma cell line Rinm5f; sequencing of a rat type II interleukin-1 receptor cDNA. Eur Cytokine Netw 5(3):319-330.
- 46. Thanhäuser A, et al. (1993) Pentoxifylline: a potent inhibitor of IL-1 and IFNgamma biosynthesis and BCG-induced cytotoxicity. Immunology 80:151-156.
- Raineri I and Senn HP. (1992) HIV-1 promotor insertion revealed by selective detection of chimeric provirus-host gene transcripts. Nucleic Acids Research 20:6261-6266.
- 48. Obata F, et al. (1993) A single universal primer for the T-cell receptor (TCR) variable genes enables enzymatic amplification and direct sequencing of TCR cDNA of various T-cell clones. Human Immunology 36:163-167.
- 49. Spurkland A. (1992) Magnetic isolation of mRNA for in vitro amplification. Trends in Genetics 8(7):225-226.
- Murtagh JJ, et al. (1994) Alternative splicing of the guanine nucleotidebinding regulatory protein Goa generates four distinct mRNAs. Nucleic Acids Res. 22(5):842-849.
- 51. Tebbe CC, et al. (1995) Direct detection of recombinant gene expression by two genetically engineered Yeasts in soil on the transcriptional and translational levels. Appl Environ Microbiology 61(12):4296-4303.
- 52. Faulkner JDB and Minton NP. (1993) Rapid small-scale isolation of mRNA from whole yeast cells. Biotechniques 14(5):718-720.
- 53. Chiodi F, et al. (1992) Human immuno-deficiency virus type 1 is present in the cerebrospinal fluid of a majority of infected individuals. J Clin Microbiol 30(7):1768-1771.
- 54. Meijer A, et al. (1995) Development of molecular methods for detection and epidemiological investigation of HIV-1, HIV-2 and HTLVI/II infections. Report 118504 001 National Inst. Publ. Health and Env. Protection Bilthoven, The Netherlands.
- 55. Albert JA, et al. (1992) Persistence of azidothymidine-resistant human immunodeficiency virus type 1 RNA genotypes in posttreatment sera. J Virol 66(9):5627-5630.

- 56. Scarlatti G, et al. (1993) Comparison of variable region 3 sequences of human immunodeficiency virus type 1 from infected children with the RNA and DNA sequences of the virus population of their mothers. Proc Natl Acad Sci USA 90:1721-1725.
- 57. Sin Y, et al. (2008) Skin bradykinin-related peptides (BRPs) and their biosynthetic precursors (kininogens): Comparisons between various taxa of Chinese and North American ranid frogs. Peptides 29:393-403.
- 58. Lu X, et al. (2007) A comparison of CFU-GM, BFU-E and endothelial progenitor cells using ex vivo expansion of selected cord blood CD133+ and CD34+ cells. Cytotherapy 9:292 - 300.
- 59. Silva J, et al. (2007) Circulating Bmi-1 mRNA as a possible prognostic factor for advanced breast cancer patients. Breast Cancer Research 9:1-9.
- 60. Jost R, et al. (2007) Magnetic quantitative reverse transcription PCR: A highthroughput method for mRNA extraction and quantitative reverse transcription PCR. BioTechniques 43:206-211.

©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

SPEC-05892

Headquarters 5791 Van Allen Way | Carlsbad, CA 92008 USA Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

