

Dynabeads™ Oligo (dT)₂₅

Catalog nos. 61002, 61005

Publication No. MAN0019067

Store at 2°C to 8°C

Rev. B

Product contents

Cat. No.	Volume
61002	2 mL
61005	5 mL

Dynabeads™ Oligo (dT)₂₅ are supplied as a suspension of approximately 5 mg/mL in PBS pH 7.4 containing 0.02% sodium azide as a preservative.

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

Product description

Dynabeads™ Oligo (dT)₂₅ are designed for the rapid isolation of highly purified, intact mRNA from eukaryotic total RNA or directly from crude extracts of cells, animal and plant tissues. The isolated mRNA can be used directly in most downstream applications in molecular biology: RT-PCR, solid-phase cDNA library construction, S1 nuclease analysis, ribonuclease protection assay, primer extension, dot and slot hybridization, *in vitro* translation experiments, RACE, subtractive hybridization, northern analysis, gene cloning, and gene expression analysis. The use of Dynabeads™ Oligo (dT)₂₅ relies on base-pairing between the poly A tail of messenger RNA and the oligo dT sequences bound to the surface of the beads. After annealing, the vial is placed on a magnet (Dynal MPC™) to concentrate the beads with their bound mRNA at the side of the tube. The supernatant containing unwanted contaminants is discarded. The protocol can be performed in 15 minutes, without the need to prepare total RNA or perform any other purification steps. The oligo dT bound to the bead surface can be used to capture the mRNA and act as a primer for reverse transcriptase during first strand cDNA synthesis. As the oligo dT is covalently bound to the Dynabeads™ surface it is possible to regenerate the Dynabeads™ Oligo (dT)₂₅ for reuse.

Required materials

- DynaMag™ Magnet (See thermofisher.com/magnets for recommendations)
- Mixing device with tilting and rotation (e.g. HulaMixer™ Sample Mixer).
- Sterile and RNase-free test tubes and pipette tips.
- Buffers/solutions (see Table 1).
- Water bath or heating block.

For tissue samples only:

- Liquid nitrogen
- Manual tissue grinder
- Syringe and needle

General guidelines

- Always keep the beads in liquid suspension, since dried out beads can lead to reduced isolation efficiency.
- The complete removal of all the buffer during washing is extremely important when working with small volumes.
- If you have cells previously isolated using other Dynabeads™ products, remove all the beads from the lysate before adding Dynabeads™ Oligo (dT)₂₅.
- We recommend the bead-mRNA complex to be used immediately for RT-PCR. If storage is necessary, elute the mRNA from the beads and freeze.
- RNases are very stable, active enzymes and generally require no cofactors to function. RNase inhibitors may be added to the protocol at any step, although this is normally not necessary. If storage of the eluted mRNA is required, addition of an RNase inhibitor at the elution step is recommended.
- Minimize the activity of RNases by creating a ribonuclease-free environment.
- Wear disposable gloves at all times during the procedure. Change gloves frequently when working with RNA (particularly if they come into contact with "dirty" glassware and surfaces).

Recommended buffers/solutions

All common buffers for mRNA purification and isolation can be used with Dynabeads™ Oligo (dT)₂₅. To take full advantage of the unique properties of the beads, the buffers described in the following sections are recommended. All buffers should be brought to room temperature prior to use, apart from the 10 mM Tris-HCl which should be kept on ice or at 2°C to 8°C.

Table 1: Recommended buffers and solutions

Binding Buffer	20 mM Tris-HCl, pH 7.5, 1.0 M LiCl, 2 mM EDTA.
Lysis/Binding Buffer	100 mM Tris-HCl, pH 7.5, 500 mM LiCl, 10 mM EDTA, 1% LiDS, 5 mM dithiothreitol (DTT). If any precipitation is observed, warm the buffer to room temperature and shake until all the components are fully resuspended.
Washing Buffer A	10 mM Tris-HCl, pH 7.5, 0.15 M LiCl, 1 mM EDTA, 0.1% LiDS.
Washing Buffer B	10 mM Tris-HCl, pH 7.5, 0.15 M LiCl, 1 mM EDTA

Note: All reagents used should be analytical grade and RNase-free.

Preparation of RNase-free solutions and materials

Water and salt solutions used in RNA preparation should be RNase-free. Wherever possible, the solutions should be treated with 0.1% diethylpyrocarbonate (DEPC) for at least 1 hour at 37°C and then heated to 100°C for 15 minutes or autoclaved for 15 minutes to remove any traces of DEPC. Tris Buffers cannot be DEPC-treated, as Tris inactivates DEPC. Solutions should be DEPC-treated and autoclaved before adding Tris. After addition of Tris, the solution should be autoclaved again. DEPC is a suspected carcinogen and should be handled with care. Sterile, disposable plasticware is essentially free of RNases and can be used for the preparation and storage of RNA without pre-treatment. General laboratory plasticware should be rinsed with chloroform.

Preparation of mRNA for downstream applications

For northern analysis, the mRNA can be eluted directly into a loading buffer containing formamide and loaded directly onto the gel. If the mRNA is to be used in downstream enzymatic applications (cDNA synthesis, *in vitro* translations experiments, RT-PCR), detergents should be omitted in the final washing steps and the elution step. Enzymatic downstream applications are not inhibited by the presence of the beads. It is possible to construct solid-phase cDNA libraries specific for a particular cell type or tissue directly on the bead-surface. The covalently linked oligo dT sequence is used to capture the mRNA and as a primer for the reverse transcriptase to synthesize the first strand cDNA. This results in a covalently linked first-strand cDNA library.

Protocol

Wash Dynabeads™ magnetic beads

- Resuspend the beads in the vial (i.e. vortex for >30 sec, or tilt and rotate for 5 min).
- Transfer the desired volume of beads to a tube.
- Add the same volume of Binding Buffer, or at least 1 mL, and resuspend.
- Place the tube in a magnet for 1 min and discard the supernatant.
- Remove the tube from the magnet and resuspend the washed beads in the same volume of Binding Buffer as the initial volume of beads taken from the vial (step 2).

Prepare lysate (animal and plant tissue)

This protocol is recommended for a sample size of 20–50 mg of solid animal tissue or 100 mg plant tissue, but can be scaled up or down to suit specific sample size requirements.

- Prepare aliquots of frozen animal or plant tissue in the amount specified for the tissue type. Do not exceed the recommended amount, or mRNA yield and purity can be reduced.
- Grind frozen tissue in liquid nitrogen. Ensure tissue remains frozen at all times to avoid RNA degradation.

- 3. Transfer the frozen powder to a homogenizer containing 1 mL Lysis/Binding Buffer and homogenize for 1–2 min until the tissue has completely lysed. Rapid lysis in Lysis/Binding Buffer prevents degradation of mRNA.
- 4. (Optional) If the raw extract is noticeably viscous, reduce the viscosity by shearing the DNA. Pass the lysate through a 21 gauge needle with a 1–2 mL syringe three times. Repeated shearing can cause the lysate to foam, but should not affect the mRNA yield. The foam can be reduced by centrifugation for 30 sec.
- 5. Centrifuge the lysate for 30–60 sec in a microcentrifuge to remove debris. The lysate is now ready for mRNA isolation (see “Isolation of mRNA from crude lysate”) or can be frozen and stored at –80°C for later use.

Prepare lysate (cultured cells and cell suspensions)

This protocol is recommended for 1–4 × 10⁶ cells, but can be scaled up or down to suit specific sample size requirements.

- 1. Wash the cell suspension in phosphate-buffered saline (PBS) and centrifuge to obtain a cell pellet. The cell pellet can be used immediately, or frozen in liquid nitrogen and stored at -80°C for later use. A stored cell pellet should be used directly from frozen.
- 2. Add 1.0 mL Lysis/Binding Buffer to the cell pellet (1–4 × 10⁶ cells). Pipet up and down a couple of times to ensure complete lysis. The release of DNA during lysis results in a viscous solution which confirms complete lysis.
- 3. Reduce the viscosity by shearing the DNA. Pass the lysate through a 21 gauge needle with a 1–2 mL syringe three times. Repeated shearing can cause the lysate to foam, but should not affect the mRNA yield. The foam can be reduced by centrifugation for 30 sec.
- 4. The lysate is now ready for mRNA isolation (see "Isolation of mRNA from crude lysate") or can be frozen and stored at –80°C for later use.

Isolate mRNA from crude lysate

- 1. Remove the solution from the washed Dynabeads® Oligo (dT)₂₅ (see "Wash Dynabeads®") and add the lysate (from section "Prepare lysate from solid animal and plant tissues" or "Prepare lysate from cultural cells and cell suspensions").
- 2. Mix beads and lysate. Allow binding by rotating on a mixer for 3–5 min at room temperature, increase the annealing time if the solution is viscous. During this step the mRNA anneals to the oligo dT sequence.
- 3. Place the vial on the magnet for 2 min and remove the supernatant.
- 4. Wash the beads twice in room temperature using the magnet: Wash once with 1 mL Washing Buffer A, and once with 1 mL Washing Buffer B. Resuspend the beads thoroughly in the Washing Buffers to remove possible contaminants, and remove the supernatant completely between the washing steps.
- 5. Perform *one* of the following:
 - a. If the bead-bound isolated mRNA is to be used in enzymatic downstream applications (e.g. solid-phase cDNA synthesis), wash one extra time with Washing Buffer B (500 µL) followed by one wash with the enzymatic buffer used in the downstream application.
 - b. To elute mRNA from the beads, remove the Washing Buffer B and add 10–20 µL 10 mM Tris-HCl. Incubate at 75°C to 80°C for 2 min, then place the tube on the magnet and quickly transfer the supernatant containing the mRNA to a new RNase-free tube. The final yield may vary somewhat between tissues/cells depending on mRNA abundance.

Purify mRNA from total RNA

This protocol describes purification of mRNA from 75 µg of total RNA, but can be scaled up or down to suit specific sample size requirements.

- 1. Adjust the volume of the 75 µg total RNA sample to 100 µL with distilled DEPC-treated water or with 10 mM Tris-HCl pH 7.5.
- 2. Add 100 µL of Binding Buffer. If total RNA is more dilute than 75 µg/100 µL, then simply add an equal volume of Binding Buffer to the beads.
- 3. Heat to 65°C for 2 min to disrupt secondary structures. Immediately place on ice.
- 4. Add the 200 µL of total RNA to the 100 µL washed beads (see “Wash Dynabeads®” section). For every 75 µg total RNA, use 1 mg beads which are washed and resuspended in 100 µL of Binding Buffer.
- 5. Mix thoroughly and allow binding by rotating continuously on a mixer for 5 min at room temperature.
- 6. Place the tube on the magnet for 1–2 min and carefully remove all the supernatant.

- 7. Remove the tube from the magnet and add 200 µL Washing Buffer B. Mix by pipetting carefully a couple of times.
- 8. Apply to the magnet for 1 min and remove the supernatant.
- 9. Repeat steps 7–8 once.
- 10. Perform *one* of the following:
 - **mRNA does not need to be eluted off the beads**
wash one more time with the same buffer that will be used in the downstream application.
 - **mRNA to be eluted off the beads**
Remove the Washing Buffer B and add 10–20 µL 10 mM Tris-HCl. Incubate at 75°C to 80°C for 2 min, then place the tube on the magnet and quickly transfer the supernatant containing the mRNA to a new RNase-free tube.

Description of materials

Dynabeads® Oligo (dT)₂₅ are uniform superparamagnetic, monodisperse polymer particles with oligo dT sequences covalently coupled to the bead surface.

Binding Capacity

Up to 2 µg poly(A)⁺ RNA can be isolated per 200 µL (1 mg) of beads, depending on the tissue or cell type and the expression level of the mRNA. A typical mammalian cell contains about 10–30 pg of RNA of which 1–5% is mRNA. The total capacity per mL of beads is approx. 10 µg mRNA. If the same beads are reused for a total of 5 mRNA isolations (four regeneration cycles) the total capacity of 1 mL beads is up to 50 µg of mRNA.

Related products

Product	Cat. No.
DynaMag™-2 Magnet	12321D
DynaMag™-5 Magnet	12303D
HulaMixer™ Sample Mixer	15920D
Phosphate Buffered Saline	14190
Dynabeads™ mRNA DIRECT™ Kit	61011
Dynabeads™ mRNA DIRECT™ Micro Kit	61021
Dynabeads™ mRNA Purification Kit	61006

[REF] on labels is the symbol for catalog number.

Important licensing information

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Manufactured by Thermo Fisher Scientific AS, Ullernchausseen 52 PO Box 114, Smestad, Oslo, Norway N-0379 and by Thermo Fisher Scientific Baltics UAB, V.A. Graiciuno 8, LT-02241 Vilnius, Lithuania. Thermo Fisher Scientific AS and Thermo Fisher Scientific Baltics UAB comply with the Quality System Standards ISO 9001 and ISO 13485.

Limited product warranty

Life Technologies Corporation and its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.

Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288
© 2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.
DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

For support visit thermofisher.com/support or email techsupport@lifetech.com