

Fresh Tissue	Frozen Tissue	Cells Grown in a Monolayer	Suspension Cells
<p>Immediately add 1 ml of PureZOL™ reagent to up to 100 mg of freshly dissected tissue.</p> <p>Disrupt and homogenize vigorously for 30–60 sec.</p> <p>Incubate lysate at room temp. for 5 min.</p>	<p>Grind frozen tissue to a fine powder with a mortar and pestle under liquid nitrogen.</p> <p>Do not let the tissue thaw.</p> <p>Add 1 ml of PureZOL to up to 100 mg of tissue.</p> <p>Disrupt and homogenize vigorously for 30–60 sec.</p> <p>Incubate lysate at room temp. for 5 min.</p>	<p>Aspirate the culture medium from a 10<sup>2</sup> cm plate containing cells grown in a monolayer.</p> <p>Do not wash cells prior to adding PureZOL.</p> <p>Immediately add 1 ml of PureZOL directly in the culture dish and pipet up and down to lyse.</p> <p>Incubate lysate at room temp. for 5 min.</p>	<p>Transfer up to 1 x 10<sup>7</sup> cells into a tube and spin for 2 min.</p> <p>Do not wash cells prior to adding PureZOL.</p> <p>Immediately add 1 ml of PureZOL directly in the tube and pipet up and down to lyse.</p> <p>Incubate lysate at room temp. for 5 min.</p>

Continue with the following steps for all sample types:

**Add 0.2 ml of chloroform.** Cover and shake vigorously for 15 sec. Incubate for 5 min at room temp.

**Centrifuge at 12,000 x g for 15 min at 4°C.** Carefully transfer only the aqueous phase to a new 2.0 ml tube.

**Add an equal volume (approx. 600 µl) of 70% ethanol.** Mix thoroughly by pipetting up and down.

*Assemble manifold properly for isolation.*

**Transfer 700 µl of lysate.** Apply vacuum. Repeat with remaining lysate.

**Add 700 µl low-stringency wash.** Apply vacuum.

*Dilute 5 µl reconstituted\* DNase I with 75 µl DNase dilution solution.*

**Add 80 µl diluted DNase I.** Incubate 15 min at room temp. Apply vacuum.

**Add 700 µl high-stringency wash.** Apply vacuum.

**Add 700 µl low-stringency wash.** Apply vacuum. Spin-purge 2 min into a 2 ml capless tube.

*Place RNA binding column into a 1.5 ml capped tube.*

**Add 40 µl (or 30 µl)\*\* of 70°C elution solution onto the center of the membrane stack.** Incubate 1 min. Centrifuge 2 min to elute.

**Repeat the elution with another 40 µl\*\* of elution solution.**

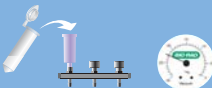
0.2 ml chloroform



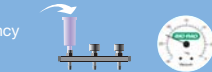
70% ethanol



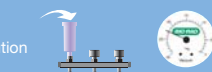
700 µl Lysate



700 µl low-stringency wash



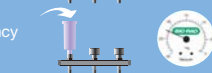
80 µl DNase I in dilution solution



700 µl high stringency wash



700 µl low stringency wash



40 µl (or 30 µl)\*\* elution solution (repeat if necessary)



\* Refer to manual for detailed protocol.

\*\* When isolating total RNA from small amounts of starting material (<10 mg of tissue or 500,000 cells), perform a single elution with 30 µl of warmed elution solution. Do not repeat the second elution.

**BIO-RAD**

Fresh Tissue	Frozen Tissue	Cells Grown in a Monolayer	Suspension Cells
<p>Immediately add 1 ml of PureZOL™ reagent to up to 100 mg of freshly dissected tissue.</p> <p>Disrupt and homogenize vigorously for 30–60 sec.</p> <p>Incubate lysate at room temp. for 5 min.</p>	<p>Grind frozen tissue to a fine powder with a mortar and pestle under liquid nitrogen.</p> <p>Do not let the tissue thaw.</p> <p>Add 1 ml of PureZOL to up to 100 mg of tissue.</p> <p>Disrupt and homogenize vigorously for 30–60 sec.</p> <p>Incubate lysate at room temp. for 5 min.</p>	<p>Aspirate the culture medium from a 10<sup>2</sup> cm plate containing cells grown in a monolayer.</p> <p>Do not wash cells prior to adding PureZOL.</p> <p>Immediately add 1 ml of PureZOL directly in the culture dish and pipet up and down to lyse.</p> <p>Incubate lysate at room temp. for 5 min.</p>	<p>Transfer up to 1 x 10<sup>7</sup> cells into a tube and spin for 2 min.</p> <p>Do not wash cells prior to adding PureZOL.</p> <p>Immediately add 1 ml of PureZOL directly in the tube and pipet up and down to lyse.</p> <p>Incubate lysate at room temp. for 5 min.</p>

Continue with the following steps for all sample types:

**Add 0.2 ml of chloroform.** Cover and shake vigorously for 15 sec. Incubate for 5 min at room temp.

**Centrifuge at 12,000 x g for 15 min at 4°C.** Carefully transfer only the aqueous phase to a new 2.0 ml tube.

**Add an equal volume (approx. 600 µl) of 70% ethanol.** Mix thoroughly by pipetting up and down.

*Insert RNA binding column into a 2 ml capped tube.*

**Transfer 700 µl of lysate.** Centrifuge 30 sec. Discard filtrate. Repeat with remaining lysate.

**Add 700 µl low-stringency wash.** Centrifuge 30 sec. Discard filtrate.

*Dilute 5 µl reconstituted\* DNase I with 75 µl DNase dilution solution.*

**Add 80 µl diluted DNase I.** Incubate 15 min at room temp. Centrifuge 30 sec. Discard filtrate.

**Add 700 µl high-stringency wash.** Centrifuge 30 sec. Discard filtrate.

**Add 700 µl low-stringency wash.** Centrifuge 1 min. Discard filtrate. Centrifuge additional 2 min.

*Place RNA binding column into a 1.5 ml capped tube.*

**Add 40 µl (or 30 µl)\*\* of 70°C elution solution onto the center of the membrane stack.** Incubate 1 min. Centrifuge 2 min to elute.

**Repeat the elution with another 40 µl\*\* of elution solution.**

0.2 ml chloroform



70% ethanol



700 µl Lysate



700 µl low-stringency wash



80 µl DNase I in dilution solution



700 µl high-stringency wash



700 µl low-stringency wash



40 µl (or 30 µl)\*\* elution solution (repeat if necessary)



\* Refer to manual for detailed protocol.

\*\* When isolating total RNA from small amounts of starting material (<10 mg of tissue or 500,000 cells), perform a single elution with 30 µl of warmed elution solution. Do not repeat the second elution.