









| Cultured cells   |  | Bacterial cells   | Yeast cells  |
|--|--|---|--|
| <p><b>Adherent</b><br/>Rinse vessel with PBS, aspirate.<br/>Lyse in vessel if # of cells &lt;math&gt;2 \times 10^6&lt;/math&gt;.</p> | <p><b>Nonadherent</b><br/>Rinse with PBS.<br/>Transfer up to <math>2 \times 10^6</math> cells, centrifuge 2 min.<br/>Decant supernatant.</p> | <p>Transfer up to <math>2.4 \times 10^9</math> cells into a capped 2 ml tube. Centrifuge at maximum speed 1 min.<br/>Decant supernatant.<br/>Add 100 <math>\mu</math>l of 500 <math>\mu</math>g/ml lysozyme. Pipet up and down.<br/>Incubate at room temp. for 5 min.</p> | <p>Transfer up to <math>3 \times 10^7</math> cells into a capped 2 ml tube. Centrifuge at maximum speed 1 min.<br/>Decant supernatant.<br/>Add 1 ml of 50 U/ml lyticase in lyticase dilution buffer. Pipet up and down.<br/>Incubate at room temp. for 10 min. Centrifuge at 5,000 rpm for 5 min. Discard supernatant.</p> |
| <p>Add 350 <math>\mu</math>l lysis solution. Pipet up and down 12x.</p>  | <p>Add 350 <math>\mu</math>l 70% EtOH. Pipet up and down.</p>  | <p>Add 350 <math>\mu</math>l lysis solution. Pipet up and down 12x.</p>   | <p>Add 350 <math>\mu</math>l 70% EtOH. Pipet up and down.</p>  |
|   |   |    |   |
|   |   |    |   |

Continue with the following steps for all sample types:



**Assemble manifold properly for isolation.**

**Transfer lysate to RNA binding column.**

Apply vacuum.

Homogenized lysate



**Add 700  $\mu$ l low stringency wash.**

Apply vacuum.

700  $\mu$ l low stringency wash



**Dilute 5  $\mu$ l reconstituted\* DNase I with 75  $\mu$ l DNase dilution solution.**

**Add 80  $\mu$ l diluted DNase I.**

Incubate 15 min at room temp. Apply vacuum.

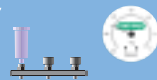
80  $\mu$ l DNase I in dilution solution



**Add 700  $\mu$ l high stringency wash.**

Apply vacuum.

700  $\mu$ l high stringency wash



**Add 700  $\mu$ l low stringency wash.**

Apply vacuum. Spin-purge 2 min into a capless 2 ml tube.

700  $\mu$ l low stringency wash



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

**Add 80  $\mu$ l 70°C elution solution onto membrane stack.**

Incubate 1 min. Centrifuge 2 min to elute.

80  $\mu$ l elution solution



\* Refer to manual for detailed protocol.

## Animal tissue

Cut tissue into small pieces (<5 mm).  
Grind into fine powder under liquid nitrogen.  
Do not let tissue thaw.

Transfer up to 20 mg (hard tissue) or  
40 mg (soft tissue) to a capped 2 ml tube.



## Plant tissue

Cut tissue into small pieces (<5 mm).  
Grind into fine powder under liquid nitrogen.  
Do not let tissue thaw.

Transfer up to 60 mg to  
a capped 2 ml tube.

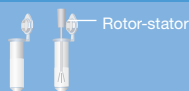


Continue with the following steps for all sample types:

### Add 700 µl lysis solution.

Disrupt vigorously with rotor-stator for 30–60 sec.

700 µl  
lysis solution



### Centrifuge lysate at maximum speed for 3 min.

Transfer supernatant to a new 2 ml capped tube.



### Add 700 µl EtOH (use 60% for animal tissue, 70% for plant tissue) to supernatant.

Homogenize with rotor-stator 30 sec.

700 µl  
60% EtOH  
or  
70% EtOH



### Assemble manifold properly for isolation.

**Transfer lysate.**  
Apply vacuum.

Homogenized  
lysate



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

700 µl  
low stringency  
wash



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

### Add 80 µl diluted DNase I.

Incubate at room temp. 25 min for animal tissue, 15 min for plant tissue. Apply vacuum.

80 µl  
DNase I in  
dilution  
solution



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

700 µl  
high stringency  
wash



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

700 µl  
low stringency  
wash



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



**Add 80 µl 70°C elution solution onto  
membrane stack.**  
Incubate 1 min. Centrifuge 2 min to elute.

80 µl  
elution  
solution



\* Refer to manual for detailed protocol.

