







Cultured cells		Bacterial cells	Yeast cells
<p>Adherent Rinse vessel with PBS, aspirate. Lyse in vessel if # of cells <2 x 10⁶.</p>	<p>Nonadherent Rinse with PBS. Transfer up to 2 x 10⁶ cells, centrifuge 2 min. Decant supernatant.</p>	<p>Transfer up to 2.4 x 10⁹ cells into a capped 2 ml tube. Centrifuge at maximum speed 1 min. Decant supernatant. Add 100 µl of 500 µg/ml lysozyme. Pipet up and down. Incubate at room temp. for 5 min.</p>	<p>Transfer up to 3 x 10⁷ cells into a capped 2 ml tube. Centrifuge at maximum speed 1 min. Decant supernatant. Add 1 ml of 50 U/ml lyticase in lyticase dilution buffer. Pipet up and down. Incubate at room temp. for 10 min. Centrifuge at 5,000 rpm for 5 min. Discard supernatant.</p>
<p>Add 350 µl lysis solution. Pipet up and down 12x.</p> 		<p>Add 350 µl lysis solution. Pipet up and down 12x.</p> 	<p>Add 350 µl lysis solution. Pipet up and down 12x.</p> 
<p>Add 350 µl 70% EtOH. Pipet up and down.</p> 		<p>Add 250 µl 70% isopropyl alcohol. Pipet up and down.</p> 	<p>Add 350 µl 70% EtOH. Pipet up and down.</p> 

Continue with the following steps for all sample types:

Insert RNA binding column into a 2 ml capless tube.

Transfer lysate.

Centrifuge 30 sec. Discard filtrate.

Homogenized lysate



Add 700 µl low stringency wash.

Centrifuge 30 sec. Discard filtrate.

700 µl low stringency wash



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.

80 µl DNase I in dilution solution



Add 700 µl high stringency wash.

Centrifuge 30 sec. Discard filtrate.

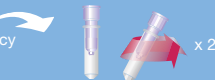
700 µl high stringency wash



Add 700 µl low stringency wash.

Centrifuge 1 min. Discard filtrate.
Centrifuge additional 2 min.

700 µl low stringency wash

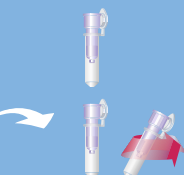


Place RNA binding column into a 1.5 ml capped tube.

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.

Aurum Total RNA Mini Kit: Cat. #732-6820

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.

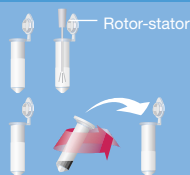


Continue with the following steps for all sample types:

Add 700 µl lysis solution.

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

700 µl
lysis solution



Centrifuge lysate at maximum speed 3 min.

Transfer supernatant to a new 2 ml capped tube.

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

Homogenize with rotor-stator 30 sec.

700 µl
60% EtOH
or
70% EtOH



Insert RNA binding column into a 2 ml capless tube.

Transfer lysate, centrifuge 60 sec.

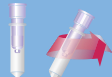
Discard filtrate. Repeat if necessary.



Add 700 µl low stringency wash.

Centrifuge 30 sec. Discard filtrate.

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. Centrifuge column 30 sec.
Discard filtrate.

80 µl
DNase in
dilution
solution



Add 700 µl high stringency wash.

Centrifuge 30 sec. Discard filtrate.

700 µl
high stringency
wash



Add 700 µl low stringency wash.

Centrifuge 30 sec. Discard filtrate.

Centrifuge additional 1 min.

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.