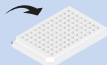


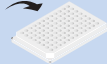


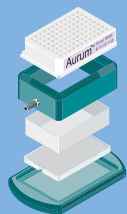


Cultured cells		Bacterial cells	Yeast cells
<b>Adherent</b> Rinse vessel with PBS, aspirate.  Lyse in vessel if # of cells $< 1 \times 10^6$ .	<b>Nonadherent</b> Rinse with PBS.  Transfer up to $1 \times 10^6$ cells per microplate well.  Centrifuge plate at $300 \times g$ for 5 min.  Aspirate supernatant.	Transfer up to $8 \times 10^8$ cells into each well of a grow block.  Centrifuge at $1,500 \times g$ for 10 min. Decant supernatant.  Add $100 \mu\text{l}$ of $500 \mu\text{g/ml}$ lysozyme. Pipet up and down.  Incubate at room temp. for 5 min.	Transfer up to $2 \times 10^7$ cells into each well of a grow block.  Centrifuge at $1,500 \times g$ for 10 min. Decant supernatant.  Add $1 \text{ ml}$ of $50 \text{ U/ml}$ lyticase in lyticase dilution buffer. Pipet up and down.  Incubate at room temp. for 10 min. Centrifuge at $1,500 \times g$ for 5 min. Discard supernatant.
Add $150 \mu\text{l}$ lysis solution. Pipet up and down. 		Add $350 \mu\text{l}$ lysis solution. Pipet up and down. 	Add $350 \mu\text{l}$ lysis solution. Pipet up and down. 
Add $150 \mu\text{l}$ 70% EtOH. Pipet up and down. 		Add $250 \mu\text{l}$ 70% isopropyl alcohol. Pipet up and down. 	Add $350 \mu\text{l}$ 70% EtOH. Pipet up and down. 

Continue with the following steps for all sample types:



**Assemble manifold properly for isolation.**

**Transfer lysate to wells of RNA binding plate.**  
Apply vacuum.

**Add  $700 \mu\text{l}$  low stringency wash to each well.**  
Apply vacuum.

**For each well, dilute  $2.5 \mu\text{l}$  reconstituted\* DNase I with  $77.5 \mu\text{l}$  DNase I dilution solution.**

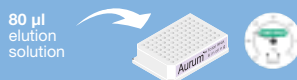
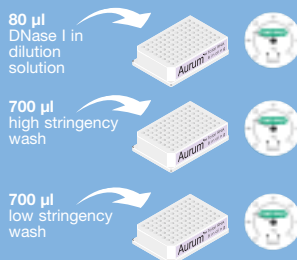
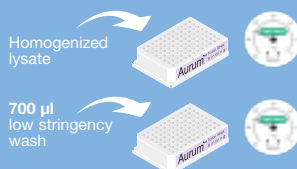
**Add  $80 \mu\text{l}$  of diluted DNase I to each well.**  
Incubate 10 min at room temp. Apply vacuum.

**Add  $700 \mu\text{l}$  high stringency wash to each well.**  
Apply vacuum.

**Add  $700 \mu\text{l}$  low stringency wash to each well.**  
Apply vacuum.

**Assemble manifold properly for elution.**

**Add  $80 \mu\text{l}$   $70^\circ\text{C}$  elution solution per well onto membrane stack.**  
Incubate 1 min. Apply vacuum gradually to  $-20$  to  $-23'' \text{ Hg}$ . Continue to apply vacuum for 5 min to elute.



\* Refer to manual for detailed protocol.

Aurum Total RNA 96 Kit: Cat. #732-6800

**BIO-RAD**