

# DNA/RNA Purification

## Selection Guide

### RNA Purification Systems

Specifications	RMag <sup>MS</sup> Total RNA	SV Total RNA	PureYield <sup>TM</sup> RNA Multiprep	SV 96 Total RNA	MagneS <sup>SV</sup> Total RNA	PolyATtract <sup>®</sup> System 1000	PolyATtract <sup>®</sup> mRNA
	Total RNA				mRNA		
	Single-Column		High-Throughput		Single		
Yields Up To (µg)	1.3–8.3 µg/mg tissue	0.27–5.3 µg/mg tissue	1mg	0.33–0.45µg/ 10 <sup>6</sup> cells	2µg/10 <sup>5</sup> cells	70 to 20ng/mg tissue	1-2% of Total RNA
Approx. Time (in minutes)	180	60	90	30	40	45	45
Technology	Sol	Mem	Mem	Mem	Mag	Mag	Mag
Manual/Automated	M	M	M	M/A	M/A	M	M
<b>Sample Type</b>							
Animal Tissues	●	●	●	●	●	●	
Cells	●	●	●	●	●	●	
Whole Blood	●	●	○		●		
Bacteria	●	●	●				
Yeast	●	●					
Plants	●	●	○				
Total RNA							●
RNA Clean-Up and Concentration		●	●				
Cat. #	Z5110	Z3100	Z3740 Z3741	Z3500 Z3505	Z3351	Z5420	Z5200

Mem – Membrane Mag – Magnetic Res – Resin Sol – Solution

● - Recommended ○ - Can be used, but may not be optimal



Promega Corporation • 2800 Woods Hollow Road • Madison, WI 53711-5399 USA • Telephone 608-274-4330 • Fax 608-277-2601  
[www.promega.com](http://www.promega.com)

©2007 Promega Corporation. All Rights Reserved.  
 Prices and specifications subject to change without prior notice.

Printed in USA, Rev 10/07

