

# DNA/RNA Purification

## Selection Guide

### Genomic DNA Purification Systems

	Wizard® Genomic DNA	Wizard® SV Genomic DNA	ReadyAmp® Genomic DNA	Wizard® SV 96 Genomic DNA	MagneSyl® Genomic Fixed Tissue	Wizard® Magnetic 96 DNA Plant	Wizard® Magnetic Food Systems	Maxwell® 16 Systems	MagneSyl® Blood Max Yield	MagneSyl® Genomic Volume Large	MagneSyl® ONE Fixed Yield	MagneSyl® Genomic KF System
	Single-Column			High-Throughput								
Specifications												
Yields Up To (µg)	Varies <sup>(1)</sup>	30	N/A	30	400ng	Varies <sup>(2)</sup>	Varies <sup>(1)</sup>	30	9	300µg/ 10ml Blood	1	6
Approx. Time (in minutes)	60	20	20	30	45	30	1 hour	40	90	8hr/96 <sup>(3)</sup>	90	25
Technology	Sol	Mem	Res	Mem	Mag	Mag	Mag	Mag	Mag	Mag	Mag	Mag
Manual/Automated	M	M	M	M/A	M	M/A	M	A	A	A	A	A
Sample Type												
Animal Blood	●		●					●	●	●	●	●
Animal Buccal Swabs		○							○		○	○
Animal Dried Blood Spots			●						○		○	○
Animal Ear Punches/Rodent Tail	●	●						●	○		○	○
Animal Stool	○						○					○
Animal Tissue	●	●		●				●	○		○	○
Bacteria (Gram-negative and -positive)	●							●				
Cells	●	●		●				●	○		○	○
Fixed Tissue (formalin, paraffin)					●							
Food							●					
Fungi	●											
Insects	●											
Plant Tissue	●	●		○		●						
Yeast	●											
Cat. #	A1120 A1125	A2360 A2361	A7710	A2370 A2371	MD1490	FF3760 FF3761	FF3750 FF3751	AS1000 AS1010 AS1020 AS1030	MD1360	A4080 A4082 A4085	MD1370	MD1460

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

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

<sup>(1)</sup> Depends on amount and sample type

<sup>(2)</sup> 100ng–2µg range

<sup>(3)</sup> Automated on Tecan Freedom EVO™ Liquid Handler



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



Promega