

DNA/RNA Purification

Selection Guide

Plasmid DNA Purification Systems

	Wizard® Plus SV Minipreps	Wizard® Plus Minipreps	Wizard® SV 96/9600	Wizard® Magnetic Bead Plasmid	Wizard® Magnetic Bead 96/9600	PureYield™ Midipreps	PureYield™ Maxipreps	Wizard® Plus Midipreps	Wizard® Plus Maxipreps	Wizard® Plus Megapreps
	Small-Scale Plasmid Systems (<20µg)					Large-Scale Plasmid Systems				
	Single-Column		High-Throughput			Single-Column				
Specifications										
Yields Up To (µg)	20	10	5	9	20	200	1mg	200	1mg	>1mg
Starting Culture Volume (ml)	1-10	1-10	1-1.5	0.5-1.5	0.5-1.5	50-250	250-500	10-100	100-500	600-1L
Approx. Time (in minutes)	45	45	40-60	45	45	30	75	90	240	240
Technology	Mem	Res	Mem	Mag	Mag	Mem	Mem	Res	Res	Res
Manual/Automated	M	M	M/A	A	A	M	M	M	M	M
Applications										
Transfection					●	●	●			
In vitro Transcription	●	●	●	●	●	●	●	●	●	●
Cloning	●	●	●	●	●	●	●	●	●	●
Probe Generation	●	●	●	●	●	●	●	●	●	●
Sequencing	●	●	●	●	●	●	●	●	●	●
PCR ^(a)	●	●	●	●	●	●	●	●	●	●
Restriction Digestion	●	●	●	●	●	●	●	●	●	●
Transformation	●	●	●	●	●	●	●	●	●	●
Cat. #	A1330 A1460	A7100 A7500	A2250 A2255 A2258	A1630 A1631	A2380 A2381	A2492 A2495	A2392	A7640	A7270	A7300

Mem – Membrane Mag – Magnetic Res – Resin
● – Recommended

^(a)The PCR process is covered by patents issued and applicable in certain countries*. Promega does not encourage or support the unauthorized or unlicensed use of the PCR process.

*In Europe, effective March 28, 2006, European Pat. Nos. 201,184 and 200,362, will expire. In the U.S., the patents covering the foundational PCR process expired on March 29, 2005.



Promega Corporation • 2800 Woods Hollow Road • Madison, WI 53711-5399 USA • Telephone 608-274-4330 • Fax 608-277-2601
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