

DNA/RNA Purification

Selection Guide

DNA Fragment Purification Systems

	Wizard® SV Gel and PCR Clean-Up	Wizard® PCR Preps	Wizard® DNA Clean-Up	Wizard® SV 96 PCR Clean-Up	Wizard® MagneSyl® PCR Clean-Up	Wizard® MagneSyl® Sequencing Reaction Clean-Up
	Single-Reaction			High-Throughput		
Specifications						
Recovery Up To	95%	96%	80%	90%	85%	Phred 20 out to at least 650 bases
Approx. Time (in minutes)	15	15	15	20	30	35
Technology	Mem	Res	Res	Mem	Mag	Mag
Manual/Automated	M	M	M	M/A	A	A
Applications						
PCR Clean-Up	●	●		●	●	
Gel Extraction	●	●		●		
Clean-up from Enzymatic Reactions	●		●			
Nucleotide Removal	●	●	●	●	●	●
Cat. #	A9281 A9282	A7170 A2180	A7280	A9341 A9342	A1930 A1931	A1831 A1832

Mem – Membrane Mag – Magnetic Res – Resin

● – Recommended



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