

DNA Analysis

Notebook

Table of Contents

Genomic DNA Purification	2	PCR Clean-Up	28
Overview		Overview	
Wizard® Genomic DNA Purification Kit		Wizard® SV Gel and PCR Clean-Up System	
Wizard® SV and SV 96 Genomic DNA Purification Systems		Wizard® SV 96 PCR Clean-Up System	
MagneSil® Blood Genomic, Max Yield System		Wizard® MagneSil® PCR Clean-Up System	
MagneSil® ONE, Fixed Yield Blood Genomic System		Wizard® PCR Preps DNA Purification System	
MagneSil® Genomic, Fixed Tissue System		Ordering Information	
MagneSil® KF, Genomic System			
Wizard® Magnetic 96 DNA Plant System			
Ordering Information			
Amplifying DNA	13	Cloning PCR DNA	34
Overview		Overview	
Optimization of PCR		Basic Subcloning	
Routine PCR		Direct Mammalian Expression	
Proofreading Polymerases		Direct Bacterial Expression	
Hot Start Methodology		PCR Cloning Techniques	
dNTPs		Ordering Information	
Ordering Information			
		DNA Analysis Tools	43
		Benchmark DNA Markers	
		Ladders and Digest Markers	