

# Introducing the Wizard® SV and SV 96 Genomic DNA Purification Systems

---

## ABSTRACT

*This article introduces the new Wizard® SV and SV 96 Genomic DNA Purification Systems. These systems can be used to purify genomic DNA from mouse tail clippings, tissue, and tissue culture cells. We have developed the Wizard® SV 96 Genomic DNA Purification System to meet the need for high-throughput genomic DNA purification. This system can be used either in a manual format or automated on a liquid handling system such as the Beckman Biomek® 2000 workstation. The Wizard® SV Genomic DNA Purification System is an easy and versatile spin or vacuum ("SV") format that provides a fast and simple way to isolate purified and intact genomic DNA in as little as 20 minutes. Genomic DNA can be purified from a variety of sample types with no detectable cross-contamination during the isolation procedure. Both systems provide similar yields of high-quality genomic DNA.*

*Promega Notes 81, 9–13.*

**Terri Grunst and Tracy Worzella**

*Promega Corporation*

Publication Date: 2002



## Download Article

(239 KB)

---