

Introducing the Maxwell® 16 Instrument: A Simple, Robust and Flexible Tool for DNA...

Introducing the Maxwell® 16 Instrument: A Simple, Robust and Flexible Tool for DNA Purification

Article

Comments & Ratings

ABSTRACT

The Maxwell® 16 Purification System combines compact instrumentation, prefilled Reagent Cartridges and optimized automated methods to maximize performance and flexibility while minimizing hands-on time required for DNA purification. The instrument is designed to process up to 16 samples in a single run of about 30 minutes. The purified DNA is ready for downstream analysis without sample preprocessing, centrifugation, precipitation or rehydration of a DNA pellet. In this article, we discuss the development of the Maxwell® 16 System and demonstrate its use for purification of genomic DNA.

Promega Notes 92, 20-23.

Dan Kephart, Steve Krueger, Terri Grunst and Hemanth Shenoi Promega Corporation

Publication Date: 2006



Download Article

(239 KB)



It appears that you have Javascript disabled. Our website requires Javascript to function correctly. For the best browsing experience, please enable Javascript.