

## Maxwell® 16 Buffy Coat Genomic DNA Purification Application

## **ABSTRACT**

The Maxwell® 16 Integrated System combines compact instrumentation, prefilled reagent cartridges and optimized automated methods to maximize performance and flexibility while minimizing hands-on time required for DNA extraction. The Maxwell® 16 Instrument can process up to 16 samples in about 30 minutes. The purified DNA is ready for downstream analysis without sample preprocessing, precipitation or rehydration of a DNA pellet. In this article, we describe extraction of DNA from buffy coat samples, the white blood cell fraction of whole blood, using the Maxwell® 16 Blood DNA Purification Kit.

Promega Notes 97, 6-8.

Hemanth Shenoi, Dan Kephart, Michael Bjerke, Sarah Shultz and Cris Cowan Promega Corporation

Publication Date: 2007



**Download Article** 

(233 KB)