

MAXWELL® 16

The Maxwell® 16 Applications Database

The Maxwell® 16 System offers automated walkaway DNA purification. The simple design of the Maxwell® 16 System and robust chemistry in the pre-filled reagent cartridges allows processing of a wide variety of sample types using the DNA IQ™ Reference Sample Kit for Maxwell® 16^(a-c) (Cat.# AS1040) or DNA IQ™ Casework Sample Kit for Maxwell® 16^(a) (Cat.# AS1210). Here we introduce the Maxwell® 16 Applications Database to facilitate sharing of Maxwell® 16-related knowledge. The database can be accessed at: www.promega.com/maxwell16/applications/ and contains information about sample types tested by both Promega scientists and Maxwell® 16 users throughout the world.

Search by keyword or browse through folders to find information about samples processed, typical yields achieved and downstream applications tested. The Forensic Applications folder contains information about DNA isolation from blood stains on a variety of solid supports, swabs, whole blood, cigarette butts and other sample types encountered in a forensic environment (Figure 1). Additional sample types are listed in the *DNA IQ™ Casework Sample Kit for Maxwell® 16 Technical Bulletin #TB354* and *DNA IQ™ Reference Sample Kit for Maxwell® 16 Technical Bulletin #TB347*.

Have you used the Maxwell® 16 Instrument to process a sample type that might be of interest to others? We encourage you to submit your Maxwell® 16 data today at: www.promega.com/maxwell16/applications/submission.asp

Visit the Maxwell® 16 Applications Database to obtain or share information about automated DNA purification from reference and casework samples.

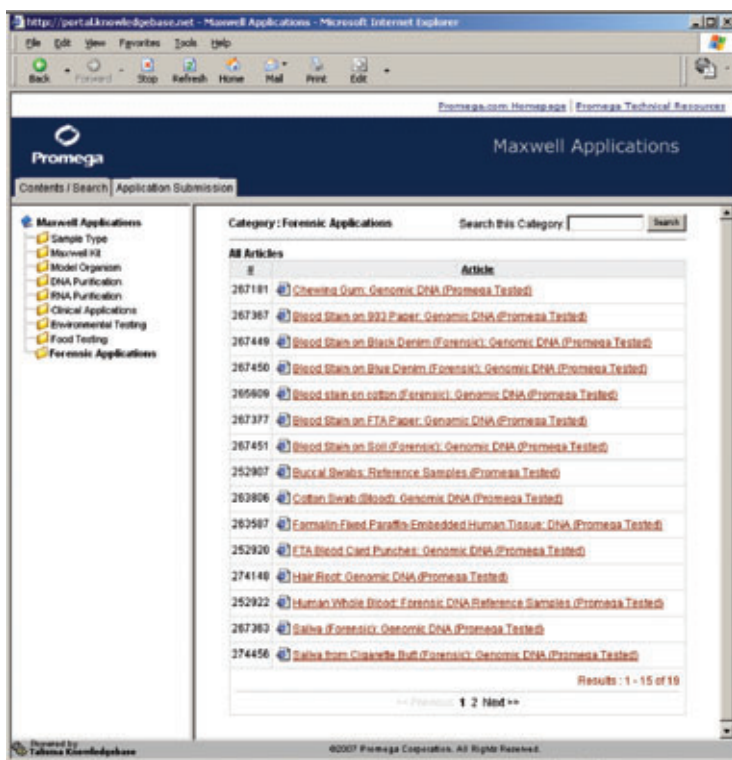


Figure 1. The Maxwell® 16 Applications Database.