

Personal Automation™ for Increased Productivity

ABSTRACT

An introduction to personal automation and how the Maxwell® 16 System provides a simple and cost-effective solution for individuals and small labs.

Promega Notes 97, 3.

Hemanth Sheno
Promega Corporation
Publication Date: 2007



Download Article

(88 KB)
