

## AccessQuick™ RT PCR System Simple Stable and Sensitive

## **ABSTRACT**

Promega's new AccessQuick $^{\text{TM}}$  RT-PCR System provides a convenient, one-step method for coupled RT-PCR. The AccessQuick $^{\text{TM}}$  Master Mix, a stable 2X solution, contains all necessary reaction components except primers, template RNA and AMV-RT, reducing setup time and pipetting steps. In testing diverse target RNAs, we found that the AccessQuick $^{\text{TM}}$  RT-PCR System provides the sensitivity and yield that researchers have today in the Access RT-PCR System.

Promega Notes 79, 12-14.

Kimberly Knoche and Mark Denhart Promega Corporation Publication Date: 2001



**Download Article** 

(183 KB)