

Plexor[®] Technology: A New Chemistry for RealTime PCR

ABSTRACT

The power of the polymerase chain reaction (PCR) is greatly enhanced by techniques that accurately, sensitively and reproducibly quantify DNA samples. These real-time or quantitative PCR (qPCR) techniques rely on the ability to detect the PCR product at each cycle during the exponential phase. Here we describe the Plexor[®] technology, a novel technique for real-time PCR that requires only two primers for sensitive and specific quantification.

Promega Notes 90, 2–4.

Susan Frackman, Steve Ekenberg, Katharine Hoffmann, Benjamin Krenke, Cynthia Sprecher and Douglas Storts

Promega Corporation

Publication Date: 2005



Download Article

(131 KB)
