

Development of a Novel Fluorescent Two-Primer Approach to Quantitative PCR

Benjamin E. Krenke, Steve Ekenberg, Susan Frackman, Katharine Hoffman, Cynthia J. Sprecher and Douglas R. Storts
Promega Corporation
Profiles in DNA 8(1), 3–5
Publication Date: 2005



Download Article

(203 KB)

Contribution of an article to *Profiles in DNA* does not constitute an endorsement of Promega products.