

High-Throughput Processing of Samples on Solid Supports Using the Slicprep™ 96 Device

Allan Tereba, Julia Krueger, Ryan Olson, Paraj Mandrekar and Bob McLaren

Promega Corporation

Profiles in DNA 8(2), 3–5

Publication Date: 2005



Download Article

(317 KB)

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