

Using Process Mapping to Improve Efficiency in a Forensic Laboratory

Julie French

Michigan State Police, Lansing, Michigan, USA

Profiles in DNA 9(2), 14–16

Publication Date: 2006



Download Article

(188 KB)

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