

# Comparison of Multiple STR Platforms and Instrumentation

---

Frank Kist, Alex Glessner, Pamela Call and Christine Tomsey

Pennsylvania State Police

*Profiles in DNA* 7(2), 3–5

Publication Date: 2004



**Download Article**

(233 KB)

---

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