

# Requirements for Complete Validation of an STR Product

---

**Thomas J. Mozer**  
Promega Corporation  
*Profiles in DNA* 4(3), 14–15  
Publication Date: 2001



## [Download Article](#)

(77 KB)

---

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