

## Y Chromosome Specific STRs

---

**Leonor Gusmão<sup>1</sup> and Angel Carracedo<sup>2</sup>**

<sup>1</sup>Instituto de Patologia e Imunologia, Molecular da Universidade do Porto, Porto, Portugal

<sup>2</sup>Institute of Legal Medicine, University of Santiago de Compostela, Santiago de Compostela, Spain

*Profiles in DNA* 6(1), 3–6

Publication Date: 2003



### Download Article

(81 KB)

---

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