

The GenePrint® PowerPlex™ 2.1 System for the FBI Selection of Thirteen CODIS Core STR...

The *GenePrint*® PowerPlex™ 2.1 System for the FBI Selection of Thirteen CODIS Core STR Loci and the Seven Standard STR Loci for ENFSI

James W. Schumm
The Bode Technology Group
Profiles in DNA 3(2), 3–7
Publication Date: 1999



Download Article

(643 KB)

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