

TWG DAM Validation of the *GenePrint® PowerPlex™ 1.1/Amelogenin and FFL Systems*

Elizabeth Amiott
Promega Corporation
Profiles in DNA 3(3), 15
Publication Date: 2000



Download Article
(368 KB)

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