DEVELOPMENT OF A NEW POWERPLEX® STR SYSTEM FOR HUMAN IDENTIFICATION

<u>Martin G. Ensenberger</u>, Patricia M. Fulmer, Robert S. McLaren, Cynthia J. Sprecher, and Douglas R. Storts

Promega Corporation, Madison, WI 53711

As human identification through STR genotyping continues to expand and mature, Promega is developing new systems to meet the evolving needs of the marketplace. Genotyping labs have increased in number and throughput; with these increases the number of profiles generated and the size of databases worldwide have grown rapidly. In turn, opportunities for data sharing and database searching are also more prevalent. Here we describe a new PowerPlex® STR system with increased discriminatory power and data sharing possibilities through the incorporation of common and informative loci used in the United States, Europe, and Asia. In addition to increased database overlap, this new system also incorporates improved chemistry to offer a robust STR solution.