

## DEVELOPMENTAL VALIDATION OF THE POWERPLEX® FUSION 6C SYSTEM

Martin Ensenberger<sup>1\*</sup>, Kristy Lenz<sup>1</sup>, Learden Matthies<sup>2</sup>, Gregory Hadinoto<sup>2</sup>, John Schienman<sup>3</sup>, Angela Przech<sup>3</sup>, Michael Morganti<sup>3</sup>, Daniel Renstrom<sup>3</sup>, Victoria Baker<sup>3</sup>, Kori Gawry<sup>4</sup>, Marlijn Hoogendoorn<sup>5</sup>, Carolyn Steffen<sup>6</sup>, Pablo Martín<sup>7</sup>, Antonio Alonso<sup>7</sup>, Hope Olson<sup>8</sup>, Cynthia Sprecher<sup>1</sup>, and Douglas Storts<sup>1</sup>

<sup>1</sup>Promega Corporation, 2800 Woods Hollow Road, Madison, WI 53711, USA

<sup>2</sup>Los Angeles County Sheriff's Department Crime Laboratory, 1800 Paseo Racho Castilla, Los Angeles, CA 90032, USA

<sup>3</sup>Connecticut Department of Emergency Services and Public Protection, Division of Scientific Services, DNA Unit, 278 Colony Street, Meriden, CT 06451, USA

<sup>4</sup>Erie County (NY) Central Police Services Forensic Laboratory, 45 Elm Street, 417E, Buffalo, NY 14203, USA

<sup>5</sup>Minnesota Bureau of Criminal Apprehension, 1430 Maryland Avenue East, St. Paul, MN 55106, USA

<sup>6</sup>National Institute of Standards and Technology, Applied Genetics Group, 100 Bureau Drive, Gaithersburg, MD 20899, USA

<sup>7</sup>National Institute of Toxicology and Forensic Sciences, Madrid Department, Biology Service, José Echegaray 4, Las Rozas 28232, Madrid, Spain

<sup>8</sup>North Dakota Office of Attorney General, Crime Laboratory Division, 2641 East Main Avenue, Bismarck, ND 58501, USA

The PowerPlex® Fusion 6C System is a 27-locus, 6-dye, multiplex that allows co-amplification and fluorescent detection of the 18 autosomal loci in the expanded CODIS core loci (CSF1PO, FGA, TH01, vWA, D1S1656, D2S1338, D2S441, D3S1358, D5S818, D7S820, D8S1179, D10S1248, D12S391, D13S317, D16S539, D18S51, D19S433 and D21S11) as well as Amelogenin and DYS391 for gender determination. The Penta D, Penta E, D22S1045, TPOX and SE33 loci are also included to increase both discrimination and overlap with databases that include profiles containing these loci. Finally, two rapidly mutating Y-STR loci, DYS570 and DYS576, are included in the multiplex. Eight Laboratories participated in the developmental validation effort following SWGDAM guidelines. Validation testing included both extracted DNA and direct amplification. Data was generated on both Applied Biosystems 3500 series and 6-dye capable 3130 series Genetic Analyzers. Results demonstrating the capabilities and limitations of the PowerPlex® Fusion 6C System are presented here and have also been submitted for NDIS review as well as for publication consideration in Forensic Science International: Genetics.