

## **DEVELOPMENTAL VALIDATION OF THE Yfiler® Plus PCR AMPLIFICATION KIT**

S. Gopinath, C. Zhong, C. Bormann Chung, V. Nguyen, M. Turnbough, J. Ge, A. Carbonaro, and J. Mulero, Thermo Fisher Scientific

Y-chromosomal markers have proven useful in solving investigations where low levels of male DNA are present in a high female DNA background. An intrinsic limitation of Y-STRs compared with autosomal STRs is a reduced power of discrimination due to a lack of recombination throughout most of the Y-chromosome. Thus, in an effort to increase the power of discrimination we have developed a new 6-dye, 27-plex Y-STR system that includes the 17 markers from the AmpF<sup>®</sup>STR Yfiler<sup>®</sup> and Yfiler Direct kits plus 10 additional highly polymorphic Y-STR markers (DYS576, DYS627, DYS460, DYS518, DYS570, DYS449, DYS481, DYF387S1a/b and DYS533). These ten new loci include 7 rapidly mutating Y-STR loci which allow for improved discrimination of related individuals.

The new multiplex is a dual application assay designed to amplify DNA from extracted casework samples and database samples from storage cards and swab lysates via direct amplification. Compared to the previous Yfiler<sup>®</sup> and Yfiler<sup>®</sup> Direct kits, the new multiplex shows improved performance in inhibited samples and admixed male and female samples at ratios >1:1000, better differentiation in male:male mixture samples in high female DNA background, and faster time to results. Additionally, no reproducible cross-reactive products were obtained on bacteria and commonly encountered animal species. The haplotype diversity and discriminatory capacity calculations for several population groups will be presented, as well as father-son studies and validation studies demonstrating improved performance with challenging samples.