

Examples of STR Population Databases For CODIS and for Casework

Bruce Budowle and Tamyra R. Moretti
FSRTC, FBI Academy, Quantico, VA 22135



The main objective for a national DNA databank is to assist investigators in the identification of perpetrators of violent crimes. For purposes of applying DNA technology to human identity testing and to make effective use of a national DNA databank, defined polymorphic genetic markers are required, and all laboratories that contribute to the database should use the same genetic loci. The short tandem repeat (STR) loci are prime candidates for typing DNA derived from forensic biological evidence and for serving as the core loci for a national DNA databank (1).

Beginning in 1996, the FBI Laboratory sponsored a community-wide science project to validate the performance of STR typing assays, to improve the quality of STR assays, to select the core STR loci to be used in the Combined DNA Index System (CODIS), and to perform population studies using the STR loci. Representatives from 21 laboratories in the United States and Canada participated in the project. The laboratories represented were the FBI, Royal Canadian Mounted Police, Armed Forces Institute of Pathology, National Institute of Standards and Technology, Alabama Department of Forensic Sciences, Arizona Department of Public Safety, California Department of Justice, Detroit Police Department, Florida Department of Law Enforcement, Illinois State Police, Metro-Dade Police Department/ Miami Children's Research Institute, Michigan State Police, Minnesota Bureau of Criminal Apprehension, North Carolina State Bureau of Investigation, Office of the Chief Medical Examiner in New York City, Orange County Sheriff's-Coroner Laboratory, Oregon State Police, Palm Beach County Sheriff's Office, Suffolk County Crime Laboratory, University of North Texas Health Science Center, and the Virginia Division of Forensic Science.

The project has been completed. As a result of this project, high quality kits that enable multiplex amplification of STR loci are commercially available, and subsequently typing procedures have been validated. Moreover, 13 core loci have been selected for use in CODIS. The loci are: CSF1PO, FGA, TH01, TPOX, vWA, D3S1358, D5S818, D7S820, D8S1179, D13S317, D16S539, D18S51, and D21S11 (1,2). The final phase of the project, which is the collection of population data, also has been completed.

Over 50 population data sets, totaling more than 10,000 individuals, comprised of African Americans, Caucasians from both the United States and Europe, Hispanics, Native Americans, and Asians have been collected and typed. These data demonstrate that the STR loci are useful for providing estimates of the frequency of a DNA profile in forensic identity testing and that a multiple locus profile is extremely rare. Furthermore, the STR loci were shown to reflect the ethnohistory of human populations, as the population data of other forensically-applied genetic markers have demonstrated. In fact, the Coefficient of Gene Differentiation based on STR loci within each population group of African Americans, Caucasians, Hispanics, and Asians is less than the NRC II Report's recommended value of 0.01 (2,3). The population data are being compiled for entry into a CODIS population data file.

To provide the reader with a general appreciation of the variation that might occur between or among major population groups, this paper provides examples of allele distributions for the 13 STR loci from each major population group relevant to the United States. Only one sample population from each major population group is displayed, because the distributions of allele frequencies for sample populations within a major population group are generally similar (except for Native Americans, where variation between Native American groups may be expected to be greater). The African American sample is from California, the Caucasian sample is from Alabama, the Hispanic sample is from Florida, the Navajo sample is from Arizona, and the Vietnamese sample was collected in California.

The distributions (in percent) of observed alleles for the 13 STR loci for five sample populations are shown in Table 1. The observed and expected homozygosities, exact test for departures from Hardy-Weinberg expectations (HWE), discrimination probability (PD), and probability of exclusion (PE) are also provided. All loci are highly polymorphic in all sample populations, with the loci D3S1358 and TH01 in Navajos having the lowest observed heterozygosities (39.6% and 48.4%, respectively), and the loci FGA in Caucasians (91.3%) and D18S51 in Hispanics (90.0%) displaying the highest observed heterozygosities. Across all sample populations, the most discriminating loci tended to be FGA, D21S11, and D18S51. The TPOX locus was the least discrimin-

ating locus in Vietnamese, Caucasians and Hispanics, and the D3S1358, D5S818, and D13S317 loci were the least discriminating loci in African Americans. These observations are similar to those reported by Budowle and Moretti (2). There was little evidence for departures from HWE in any of the populations. Based on the exact test, the loci that departed significantly from HWE are: FGA ($p=0.012$, African Americans; and $p<10^{-3}$, Navajos); TH01 ($p=0.016$, Navajos); D7S820 ($p=0.029$, Vietnamese); D8S1179 ($p=0.012$, Caucasians; and $p=0.009$, Hispanics); and D21S11 ($p=0.049$, Vietnamese; and $p=0.006$, Caucasians). After employing the Bonferroni correction for the number of loci analyzed (i.e., 13 loci per database), these observations (except, possibly that of locus FGA in Navajos) are not likely to be significant.

An inter-class correlation test analysis was performed to reveal possible correlations between alleles at any of the pair-wise comparisons of the 13 loci. For each database, there is a total of 78 pair-wise comparisons performed. The number of significant departures is at or below expected levels (5%, or 4 observations) in all groups, except Hispanics. In Hispanics, six significant departures were observed (7.7% of the pair-wise tests). However, none of the empirical levels of significance for these six observations are below the adjusted Bonferroni level. Furthermore, five of these correlations for the Hispanic sample are negative; a positive value might be attributed to the effects of substructure. Based on these observations, the data do not support any significant departure from independence between pairs of loci in any sample population. With little evidence of association between loci, the assumption of independence is valid, and a multiple-locus profile frequency can be estimated using the product rule.

Differences in allele frequencies between the loci of the major population databases displayed in Table 1 were observed at the loci and might be expected. Despite these differences in allele frequencies, a 13-locus profile frequency would be rare in all five population groups. The Navajo are the least polymorphic of the five population groups, and these data may provide insight about the

upper bound of estimates for isolated populations. The most common profile frequency derived from the 13 core STR loci is less than 1 in 950 million in the Navajo, and typically the estimates are substantially more rare. The estimates of the most common profile frequency for the other four population groups are at least two-to-three orders of magnitude smaller than that for Navajos.

In conclusion, a subset of the collected STR population data is presented. The data are similar to other relevant STR population studies and can be used for estimating the rarity of a multiple locus STR profile. Substantial population data, collected through the concerted effort of the forensic community, provide a solid base for the validity of STR typing and are available for assessing the significance of a multiple locus profile. Cooperative efforts such as the STR project are invaluable for enhancing the capabilities of forensic scientists to help resolve violent crimes.

ACKNOWLEDGEMENTS

We would like to express our thanks to those people and laboratories who contributed to and/or supported this study.

This is publication number 99-04 of the Laboratory Division of the Federal Bureau of Investigation. Names of commercial manufacturers are provided for identification only, and inclusion does not imply endorsement by the Federal Bureau of Investigation.

REFERENCES

1. Budowle B, Moretti TR, Niezgoda SJ, Brown, BL. CODIS and PCR-based short tandem repeat loci: Law enforcement tools. In: Second European Symposium on Human Identification 1998, Promega Corporation, Madison, Wisconsin (in press).
2. Budowle B, Moretti TR. Population data on the thirteen CODIS Core short tandem repeat loci in African Americans, U.S. Caucasians, Hispanics, Bahamaians, Jamaicans, and Trinidadians. J. Forens. Sci. (Submitted).
3. National Research Council II Report. The Evaluation of Forensic Evidence. National Academy Press, Washington, D.C., 1996.

Examples of STR Population Databases For CODIS and for Casework

Table 1. Observed allele distributions (as %) for 13 STR loci in five population groups (Note: these data are preliminary and may be edited slightly before being entered into CODIS)

| D3S1358 | Vietnamese (N=213) | African American (N=200) | Caucasian (N=150) | Hispanic (N=210) | Navajo (N=182) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| <12 | 0.000 | 0.250 | 0.333 | 0.238 | 0.000 |
| 12 | 0.000 | 0.500 | 0.000 | 0.000 | 0.000 |
| 13 | 0.469 | 1.250 | 0.667 | 0.952 | 1.374 |
| 14 | 3.286 | 8.500 | 15.333 | 8.095 | 3.297 |
| 15 | 30.047 | 27.500 | 23.000 | 35.238 | 73.901 |
| 16 | 34.038 | 36.250 | 25.667 | 25.238 | 13.736 |
| 17 | 24.883 | 20.500 | 22.667 | 15.714 | 6.593 |
| 18 | 5.869 | 5.000 | 10.000 | 14.286 | 0.824 |
| 19 | 1.408 | 0.250 | 2.333 | 0.238 | 0.275 |
| Homozygosity (Obs.) | 27.7% | 25.0% | 22.0% | 25.2% | 60.4% |
| Homozygosity (Exp.) | 27.1% | 25.7% | 20.2% | 23.8% | 57.0% |
| p value | 0.846 | 0.818 | 0.574 | 0.620 | 0.342 |
| exact test | 0.953 | 0.648 | 0.982 | 0.954 | 0.229 |
| PD | 0.878 | 0.889 | 0.927 | 0.905 | 0.635 |
| PE | 0.482 | 0.511 | 0.595 | 0.543 | 0.246 |

| vWA | Vietnamese (N=215) | African American (N=200) | Caucasian (N=150) | Hispanic (N=259) | Navajo (N=182) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| 13 | 0.233 | 1.250 | 0.000 | 0.386 | 0.000 |
| 14 | 26.977 | 6.250 | 10.667 | 7.336 | 2.473 |
| 15 | 1.860 | 19.250 | 9.333 | 10.232 | 1.923 |
| 16 | 14.884 | 25.250 | 21.333 | 26.641 | 42.857 |
| 17 | 22.791 | 24.500 | 28.000 | 29.923 | 31.319 |
| 18 | 23.953 | 13.250 | 20.667 | 18.533 | 15.659 |
| 19 | 7.907 | 7.250 | 8.333 | 5.405 | 5.495 |
| 20 | 1.395 | 1.500 | 1.667 | 1.544 | 0.275 |
| 21 | 0.000 | 1.000 | 0.000 | 0.000 | 0.000 |
| 22 | 0.000 | 0.500 | 0.000 | 0.000 | 0.000 |
| Homozygosity (Obs.) | 20.5% | 20.5% | 16.7% | 18.5% | 30.8% |
| Homozygosity (Exp.) | 20.9% | 18.6% | 19.1% | 21.2% | 30.8% |
| p value | 0.870 | 0.490 | 0.444 | 0.287 | 0.984 |
| exact test | 0.849 | 0.757 | 0.515 | 0.665 | 0.268 |
| PD | 0.922 | 0.936 | 0.926 | 0.920 | 0.849 |
| PE | 0.584 | 0.629 | 0.618 | 0.586 | 0.441 |

Examples of STR Population Databases For CODIS and for Casework

| FGA | Vietnamese (N=212) | African American (N=200) | Caucasian (N=150) | Hispanic (N=210) | Navajo (N=182) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| <18 | 0.943 | 0.250 | 0.000 | 0.000 | 0.000 |
| 18 | 1.887 | 0.500 | 1.333 | 0.952 | 1.374 |
| 18.2 | 0.000 | 1.500 | 0.000 | 0.000 | 0.000 |
| 19 | 8.962 | 6.750 | 4.333 | 8.333 | 18.681 |
| 19.2 | 0.000 | 0.750 | 0.000 | 0.000 | 0.000 |
| 20 | 5.896 | 4.500 | 13.667 | 12.619 | 9.066 |
| 20.2 | 0.236 | 0.250 | 0.000 | 0.000 | 0.000 |
| 21 | 15.802 | 9.500 | 17.000 | 14.524 | 12.637 |
| 21.2 | 1.415 | 0.000 | 0.000 | 0.238 | 0.000 |
| 22 | 19.340 | 17.500 | 19.000 | 14.762 | 8.516 |
| 22.2 | 0.943 | 0.250 | 1.000 | 0.476 | 0.000 |
| 23 | 12.972 | 19.750 | 14.667 | 14.048 | 5.220 |
| 23.2 | 1.887 | 0.000 | 0.333 | 0.476 | 0.000 |
| 24 | 15.094 | 16.250 | 15.667 | 14.762 | 11.264 |
| 24.2 | 1.651 | 0.000 | 0.000 | 0.000 | 0.000 |
| 25 | 6.604 | 11.250 | 11.000 | 10.714 | 16.758 |
| 25.2 | 0.943 | 0.000 | 0.000 | 0.000 | 0.000 |
| 26 | 3.538 | 4.250 | 2.000 | 4.762 | 11.264 |
| 26.2 | 0.472 | 0.000 | 0.000 | 0.000 | 0.000 |
| 27 | 1.415 | 3.500 | 0.000 | 2.619 | 4.670 |
| 28 | 0.000 | 1.750 | 0.000 | 0.476 | 0.549 |
| 30 | 0.000 | 0.000 | 0.000 | 0.238 | 0.000 |
| >30 | 0.000 | 1.500 | 0.000 | 0.000 | 0.000 |
| Homozygosity (Obs.) | 12.3% | 14.5% | 8.7% | 12.4% | 19.8% |
| Homozygosity (Exp.) | 11.9% | 12.6% | 14.2% | 12.0% | 12.3% |
| p value | 0.859 | 0.417 | 0.054 | 0.858 | 0.002 |
| exact test | 0.279 | 0.012 | 0.514 | 0.337 | 0.000 |
| PD | 0.970 | 0.965 | 0.954 | 0.969 | 0.965 |
| PE | 0.758 | 0.744 | 0.708 | 0.752 | 0.747 |

Examples of STR Population Databases For CODIS and for Casework

| TH01 | Vietnamese (N=210) | African American (N=200) | Caucasian (N=150) | Hispanic (N=240) | Navajo (N=182) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| 5 | 0.000 | 0.500 | 0.000 | 0.000 | 0.000 |
| 6 | 10.714 | 11.750 | 19.667 | 21.250 | 16.758 |
| 7 | 28.095 | 42.500 | 16.667 | 25.208 | 61.264 |
| 8 | 5.000 | 18.750 | 13.000 | 10.417 | 5.495 |
| 8.3 | 0.000 | 0.000 | 0.333 | 0.000 | 0.000 |
| 9 | 44.048 | 13.500 | 18.667 | 18.542 | 0.824 |
| 9.3 | 5.238 | 12.250 | 30.667 | 23.542 | 15.659 |
| 10 | 6.667 | 0.750 | 1.000 | 1.042 | 0.000 |
| 11 | 0.238 | 0.000 | 0.000 | 0.000 | 0.000 |
| Homozygosity (Obs.) | 32.9% | 26.0% | 20.7% | 20.0% | 51.6% |
| Homozygosity (Exp.) | 29.2% | 26.1% | 21.0% | 20.8% | 42.9% |
| p value | 0.250 | 0.973 | 0.927 | 0.766 | 0.018 |
| exact test | 0.501 | 0.313 | 0.977 | 0.390 | 0.016 |
| PD | 0.870 | 0.894 | 0.922 | 0.919 | 0.757 |
| PE | 0.478 | 0.522 | 0.582 | 0.582 | 0.339 |

| TPOX | Vietnamese (N=210) | African American (N=200) | Caucasian (N=150) | Hispanic (N=240) | Navajo (N=182) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| 6 | 0.000 | 6.250 | 0.333 | 0.417 | 0.000 |
| 7 | 0.476 | 2.500 | 0.000 | 0.208 | 0.000 |
| 8 | 56.190 | 33.250 | 54.333 | 50.625 | 34.890 |
| 9 | 9.524 | 21.000 | 13.000 | 8.333 | 0.824 |
| 10 | 6.190 | 6.000 | 4.000 | 6.250 | 1.648 |
| 11 | 25.476 | 27.000 | 24.000 | 27.708 | 37.637 |
| 12 | 2.143 | 4.000 | 4.333 | 6.458 | 25.000 |
| Homozygosity (Obs.) | 38.1% | 26.0% | 32.0% | 39.2% | 33.0% |
| Homozygosity (Exp.) | 39.3% | 23.5% | 37.1% | 34.7% | 32.4% |
| p value | 0.730 | 0.412 | 0.195 | 0.144 | 0.879 |
| exact test | 0.386 | 0.430 | 0.505 | 0.090 | 0.302 |
| PD | 0.788 | 0.908 | 0.806 | 0.829 | 0.828 |
| PE | 0.364 | 0.548 | 0.386 | 0.411 | 0.392 |

Examples of STR Population Databases For CODIS and for Casework

| CSF1PO | Vietnamese (N=210) | African American (N=200) | Caucasian (N=150) | Hispanic (N=240) | Navajo (N=182) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| 6 | 0.000 | 0.000 | 0.000 | 0.000 | 0.000 |
| 7 | 0.238 | 5.500 | 0.000 | 0.208 | 0.000 |
| 8 | 0.000 | 7.000 | 0.000 | 0.417 | 1.648 |
| 9 | 3.333 | 3.500 | 2.667 | 1.250 | 4.945 |
| 10 | 22.143 | 31.250 | 27.000 | 25.417 | 26.099 |
| 11 | 27.857 | 22.500 | 29.333 | 29.583 | 34.890 |
| 12 | 39.048 | 24.250 | 32.000 | 35.625 | 29.396 |
| 12.1 | 0.000 | 0.000 | 0.000 | 0.000 | 0.824 |
| 13 | 5.952 | 5.000 | 6.667 | 6.875 | 1.923 |
| 14 | 0.952 | 1.000 | 1.667 | 0.417 | 0.000 |
| 15 | 0.476 | 0.000 | 0.667 | 0.208 | 0.275 |
| Homozygosity (Obs.) | 24.3% | 21.0% | 26.0% | 32.1% | 28.6% |
| Homozygosity (Exp.) | 28.2% | 21.7% | 26.4% | 28.2% | 27.7% |
| p value | 0.206 | 0.813 | 0.903 | 0.187 | 0.803 |
| exact test | 0.771 | 0.999 | 0.826 | 0.675 | 0.470 |
| PD | 0.863 | 0.921 | 0.877 | 0.874 | 0.871 |
| PE | 0.471 | 0.579 | 0.490 | 0.462 | 0.471 |

| D5S818 | Vietnamese (N=212) | African American (N=200) | Caucasian (N=150) | Hispanic (N=259) | Navajo (N=182) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| 7 | 2.830 | 0.500 | 0.000 | 2.317 | 16.209 |
| 8 | 0.236 | 5.250 | 0.000 | 1.158 | 3.571 |
| 9 | 6.132 | 1.750 | 3.000 | 4.826 | 0.275 |
| 10 | 23.349 | 6.000 | 3.667 | 4.826 | 6.044 |
| 11 | 30.425 | 28.250 | 41.667 | 38.224 | 58.242 |
| 12 | 23.113 | 35.750 | 36.667 | 32.625 | 10.714 |
| 13 | 13.443 | 20.750 | 13.667 | 15.444 | 4.945 |
| 14 | 0.236 | 1.000 | 1.000 | 0.579 | 0.000 |
| 15 | 0.236 | 0.500 | 0.333 | 0.000 | 0.000 |
| >15 | 0.000 | 0.250 | 0.000 | 0.000 | 0.000 |
| Homozygosity (Obs.) | 19.8% | 26.0% | 33.3% | 30.5% | 37.9% |
| Homozygosity (Exp.) | 22.1% | 25.6% | 32.7% | 28.0% | 38.3% |
| p value | 0.416 | 0.887 | 0.865 | 0.377 | 0.922 |
| exact test | 0.096 | 0.614 | 0.673 | 0.134 | 0.837 |
| PD | 0.905 | 0.890 | 0.834 | 0.876 | 0.822 |
| PE | 0.565 | 0.515 | 0.412 | 0.483 | 0.404 |

Examples of STR Population Databases For CODIS and for Casework

| D7S820 | Vietnamese (N=212) | African American (N=200) | Caucasian (N=150) | Hispanic (N=259) | Navajo (N=182) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| 6 | 0.000 | 0.000 | 0.333 | 0.193 | 0.000 |
| 7 | 0.236 | 0.500 | 2.667 | 0.965 | 0.275 |
| 8 | 15.094 | 23.250 | 15.667 | 15.444 | 12.363 |
| 9 | 7.075 | 11.250 | 16.000 | 11.969 | 0.000 |
| 10 | 15.094 | 34.000 | 29.667 | 26.255 | 14.286 |
| 11 | 39.387 | 21.750 | 15.000 | 22.780 | 40.934 |
| 12 | 20.047 | 7.750 | 16.333 | 18.340 | 28.297 |
| 13 | 2.594 | 1.000 | 3.667 | 3.668 | 3.846 |
| 14 | 0.472 | 0.500 | 0.667 | 0.386 | 0.000 |
| Homozygosity (Obs.) | 25.9% | 26.5% | 16.0% | 19.7% | 23.6% |
| Homozygosity (Exp.) | 24.5% | 23.4% | 18.7% | 19.3% | 28.3% |
| p value | 0.621 | 0.298 | 0.401 | 0.858 | 0.163 |
| exact test | 0.029 | 0.910 | 0.675 | 0.539 | 0.029 |
| PD | 0.900 | 0.908 | 0.933 | 0.931 | 0.861 |
| PE | 0.542 | 0.547 | 0.627 | 0.614 | 0.476 |

| D18S51 | Vietnamese (N=215) | African American (N=200) | Caucasian (N=150) | Hispanic (N=210) | Navajo (N=185) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| <11 | 0.233 | 0.500 | 1.333 | 0.952 | 0.000 |
| 11 | 0.465 | 0.750 | 0.333 | 1.429 | 0.270 |
| 12 | 4.186 | 6.000 | 13.000 | 13.333 | 9.459 |
| 13 | 11.860 | 4.750 | 14.667 | 12.381 | 32.162 |
| 13.2 | 0.000 | 0.500 | 0.000 | 0.000 | 0.000 |
| 14 | 18.605 | 6.250 | 18.000 | 13.571 | 13.784 |
| 14.2 | 0.000 | 1.000 | 0.000 | 0.000 | 0.000 |
| 15 | 26.744 | 19.000 | 15.667 | 18.571 | 7.838 |
| 16 | 14.651 | 17.250 | 11.000 | 15.000 | 18.378 |
| 17 | 10.698 | 15.250 | 13.333 | 9.762 | 11.622 |
| 18 | 3.256 | 9.500 | 7.333 | 5.476 | 1.892 |
| 19 | 3.256 | 9.500 | 2.667 | 4.286 | 2.432 |
| 20 | 2.093 | 4.250 | 2.000 | 3.095 | 0.811 |
| 21 | 1.395 | 3.500 | 0.667 | 0.952 | 1.081 |
| 22 | 1.628 | 1.750 | 0.000 | 0.952 | 0.000 |
| >22 | 0.931 | 0.250 | 0.000 | 0.238 | 0.270 |
| Homozygosity (Obs.) | 19.1% | 11.0% | 16.7% | 10.0% | 20.5% |
| Homozygosity (Exp.) | 15.6% | 11.8% | 12.9% | 12.2% | 18.4% |
| p value | 0.161 | 0.717 | 0.169 | 0.325 | 0.447 |
| exact test | 0.288 | 0.363 | 0.424 | 0.957 | 0.350 |
| PD | 0.956 | 0.970 | 0.964 | 0.969 | 0.942 |
| PE | 0.690 | 0.759 | 0.732 | 0.749 | 0.643 |

Examples of STR Population Databases For CODIS and for Casework

| D8S1179 | Vietnamese (N=214) | African American (N=200) | Caucasian (N=150) | Hispanic (N=210) | Navajo (N=185) |
|---------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| <9 | 0.000 | 0.250 | 1.667 | 1.429 | 0.000 |
| 9 | 0.000 | 0.250 | 0.333 | 0.952 | 0.000 |
| 10 | 14.953 | 0.750 | 12.333 | 9.286 | 14.865 |
| 11 | 13.318 | 6.250 | 9.667 | 5.000 | 5.135 |
| 12 | 10.981 | 10.750 | 8.667 | 10.476 | 10.811 |
| 13 | 16.822 | 22.500 | 31.333 | 36.429 | 37.027 |
| 14 | 13.318 | 31.000 | 22.000 | 21.667 | 22.162 |
| 15 | 16.822 | 19.750 | 10.333 | 11.905 | 8.378 |
| 16 | 11.449 | 6.000 | 3.000 | 2.619 | 1.622 |
| 17 | 1.636 | 2.500 | 0.667 | 0.238 | 0.000 |
| 18 | 0.701 | 0.000 | 0.000 | 0.000 | 0.000 |
| Homozygosity (Obs.) | 17.8% | 22.0% | 19.3% | 27.6% | 25.4% |
| Homozygosity (Exp.) | 13.8% | 20.3% | 18.8% | 21.5% | 22.8% |
| p value | 0.092 | 0.562 | 0.863 | 0.031 | 0.395 |
| exact test | 0.052 | 0.567 | 0.012 | 0.009 | 0.682 |
| PD | 0.960 | 0.929 | 0.929 | 0.924 | 0.917 |
| PE | 0.715 | 0.602 | 0.633 | 0.595 | 0.569 |

| D16S539 | Vietnamese (N=210) | African American (N=200) | Caucasian (N=150) | Hispanic (N=240) | Navajo (N=185) |
|--------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| 5 | 0.000 | 0.250 | 0.000 | 0.000 | 0.000 |
| 8 | 0.714 | 4.000 | 0.667 | 2.292 | 0.000 |
| 9 | 22.143 | 18.250 | 11.333 | 14.583 | 16.757 |
| 10 | 16.429 | 10.500 | 5.000 | 9.583 | 16.486 |
| 11 | 26.190 | 31.000 | 31.333 | 28.125 | 14.595 |
| 12 | 23.571 | 20.750 | 31.000 | 25.417 | 35.946 |
| 13 | 9.524 | 13.000 | 19.000 | 16.250 | 14.324 |
| 14 | 1.429 | 2.250 | 1.667 | 3.542 | 1.892 |
| 15 | 0.000 | 0.000 | 0.000 | 0.208 | 0.000 |
| Homozygosity(Obs.) | 17.1% | 20.0% | 28.7% | 20.0% | 20.5% |
| Homozygosity(Exp.) | 20.8% | 20.1% | 24.4% | 20.1% | 22.5% |
| p value | 0.196 | 0.986 | 0.218 | 0.979 | 0.533 |
| exact test | 0.770 | 0.068 | 0.596 | 0.696 | 0.660 |
| PD | 0.916 | 0.923 | 0.901 | 0.929 | 0.914 |
| PE | 0.584 | 0.605 | 0.529 | 0.603 | 0.567 |

Examples of STR Population Databases For CODIS and for Casework

| D21S11 | Vietnamese (N=215) | African American (N=200) | Caucasian (N=150) | Hispanic (N=210) | Navajo (N=185) |
|--------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| <24.2 | 0.000 | 0.000 | 0.333 | 0.000 | 0.000 |
| 26 | 0.000 | 0.250 | 0.333 | 0.238 | 0.000 |
| 26.2 | 0.000 | 0.250 | 0.000 | 0.000 | 0.000 |
| 27 | 0.465 | 3.750 | 4.000 | 1.667 | 0.000 |
| 28 | 5.116 | 28.750 | 17.667 | 12.381 | 5.405 |
| 28.2 | 0.000 | 0.000 | 0.000 | 0.000 | 0.000 |
| 29 | 23.721 | 20.000 | 18.667 | 23.333 | 17.838 |
| 29.2 | 0.000 | 0.250 | 0.000 | 0.238 | 0.000 |
| 30 | 26.977 | 14.750 | 27.333 | 24.762 | 50.811 |
| 30.2 | 1.395 | 2.250 | 2.333 | 3.333 | 0.000 |
| 31 | 8.837 | 7.000 | 8.000 | 7.619 | 4.865 |
| 31.2 | 8.140 | 3.750 | 9.333 | 9.524 | 6.216 |
| 32 | 2.791 | 1.000 | 1.667 | 0.952 | 0.000 |
| 32.2 | 15.814 | 8.000 | 7.000 | 10.952 | 11.892 |
| 33 | 0.930 | 0.250 | 0.000 | 0.000 | 0.000 |
| 33.1 | 0.000 | 0.500 | 0.000 | 0.000 | 0.000 |
| 33.2 | 5.349 | 3.500 | 3.000 | 3.571 | 2.703 |
| 34 | 0.000 | 1.000 | 0.000 | 0.000 | 0.000 |
| 34.1 | 0.000 | 0.250 | 0.000 | 0.000 | 0.000 |
| 34.2 | 0.465 | 0.250 | 0.333 | 0.952 | 0.270 |
| 35 | 0.000 | 3.500 | 0.000 | 0.000 | 0.000 |
| 36 | 0.000 | 0.750 | 0.000 | 0.238 | 0.000 |
| >36 | 0.000 | 0.000 | 0.000 | 0.238 | 0.000 |
| Homozygosity(Obs.) | 20.5% | 15.0% | 18.7% | 13.3% | 32.4% |
| Homozygosity(Exp.) | 17.3% | 16.0% | 16.1% | 15.9% | 31.2% |
| P Value | 0.222 | 0.708 | 0.399 | 0.312 | 0.721 |
| Exact Test | 0.049 | 0.930 | 0.006 | 0.451 | 0.853 |
| PD | 0.946 | 0.955 | 0.944 | 0.949 | 0.875 |
| PE | 0.659 | 0.688 | 0.679 | 0.684 | 0.478 |

Examples of STR Population Databases For CODIS and for Casework

| D13S317 | Vietnamese (N=211) | African American (N=200) | Caucasian (N=150) | Hispanic (N=259) | Navajo (N=182) |
|--------------------|-----------------------|--------------------------------|----------------------|---------------------|-------------------|
| 7 | 0.000 | 0.000 | 0.000 | 0.000 | 0.000 |
| 8 | 36.256 | 4.500 | 11.667 | 11.197 | 0.824 |
| 9 | 11.374 | 3.250 | 8.000 | 11.583 | 24.725 |
| 10 | 11.611 | 2.250 | 6.667 | 7.915 | 15.110 |
| 11 | 20.142 | 27.000 | 32.000 | 31.660 | 22.253 |
| 12 | 16.825 | 40.500 | 26.667 | 22.201 | 19.780 |
| 13 | 2.133 | 15.250 | 11.000 | 10.232 | 15.934 |
| 14 | 0.948 | 7.250 | 3.667 | 5.212 | 1.374 |
| 15 | 0.711 | 0.000 | 0.333 | 0.000 | 0.000 |
| Homozygosity(Obs.) | 26.5% | 29.5% | 20.7% | 23.9% | 19.2% |
| Homozygosity(Exp.) | 22.6% | 26.7% | 20.9% | 19.3% | 19.6% |
| P Value | 0.166 | 0.374 | 0.949 | 0.061 | 0.899 |
| Exact Test | 0.940 | 0.493 | 0.389 | 0.064 | 0.799 |
| PD | 0.917 | 0.891 | 0.927 | 0.931 | 0.929 |
| PE | 0.569 | 0.508 | 0.596 | 0.624 | 0.603 |