On-Line Autosomal and Y- STR Reference Database of Argentina.

Miguel Marino, Andrea Sala. and Daniel Corach

Servicio de Huellas Digitales Genéticas y Cátedra de Genética y biología Molecular, Facultad de Farmacia y Bioquímica, Universidad de Buenos Aires, Junín 956 1113-Buenos Aires, Argentina <a href="mailto:shdq@ffyb.uba.ar">shdq@ffyb.uba.ar</a>.

Allele frequencies represent the basic tool for statistical evaluation of molecular studies based on DNA typing used in personal identification and forensic casework. In order to provide to the forensic and the molecular anthropology communities a rapid access by Internet on-line inquires a STR reference database has been organized in the website www.ffyb.uba.ar\bdg.

The database has a friendly access to the required information from each province of the entire country. On karyotype scheme it can be located each of the thirteen CODIS markers and on the Y chromosome each of the minimal haplotype STRs included initially at the <a href="www.yhrd.org">www.yhrd.org</a>. The frequencies are updated every September but previous years would be checked in order to confirm previous calculations. In addition to the database that includes over 2700 genotypes a complete list of references is included. Anonymous genotypes are included and most samples were obtained from routine casework.

At present over nine provinces are included and an additional support will be offered for aboriginal population data, mainly well defined ethnic groups typed either by autosomal and Y STRs as well as SNPs and mtDNA sequence data. Although it is complete free of charge, it requires registration for access.