

Abstracts

Two Different VNTR Loci (3' apoB and D1S80) in Yugoslav Population and their Application in Paternity Testing

Sanja Glišić, Slaven Erceg, Nataša Radovanović, Dragan Alavantic
VINCA Institute, Lab for radiob and mol genetics, p.o. box 522, 11000 Belgrade, Yugoslavia

~~~~~

We investigated polymorphism in minisatellite region caused by a variable number of tandem repeats located at the 3'end of the human apolipoprotein B gene in 696 Serbian (Belgrade area) unrelated individuals of both genders. After using the polymerase chain reaction (PCR) 17 different alleles, containing 22-54 repeats were distinguished and sixty three different 3' apoB HVR genotypes. Heterozygosity for 3' apoB HVR was 0.71. It was registered biomodal frequency distribution (picks: HVE34 and HVE44).

Techniques we applied: DNA extraction was performed by Kunkel's method for DNA isolation from whole blood and from other sources (from hair, sperm, bloodstains, biopsies from different tissues) and amplified fragment length polymorphism (AFLP) was followed obviously. HVR alleles were distinguished on 1.2% agarose gel and visualized by ethidium-bromide. The results were analyzed by use of UVP GSD8000 Gel Documentation System.

ପୁରୀ