

Population Data of the Four STR Systems: D3S1358, D8S1179, D18S51 and D21S11

Esther Mornhinweg, Christine Luckenbach, Horst Ritter



A SW- German population has been investigated for four STR systems: the three compound repeats, D3S1358 (n=500) and D18S51 (n=450) and the complex-STR D21S11 (n=500).

Separation was carried out manually on a polyacrylamide gel and automatically with the ABI 310 Genetic Analyzer.

All four systems meet Hardy-Weinberg expectations and a comparison with other populations showed no significant differences. Additionally we present the mutation rates of all four systems analyzing 65 families.