

COMPARATIVE ANALYSIS OF 10 X-STR OF A PERUVIAN POPULATION WITH 15 IBERIAN AND LATIN AMERICAN POPULATIONS

Gavilan M, Talledo M, Aiquipa L, Choque C, Almeyda ML, Arevalo J, and Montoya Y
Paternity Laboratory, BIOLINKS S.A. Javier Prado Oeste Avenue 844 Magdalena Lima 17, Peru

A study on a population of the city of Lima, which shows characteristics of ancestral Andean of Peru, was evaluated with ten X-chromosome short tandem repeats (X-STR), which included DXS8378, DXS9902, DXS7132, DXS9898, DXS6809, DXS6789, DXS7133, GATA172D05, GATA31E08, and DXS7423. Each of the markers was evaluated on 300 unrelated individuals (150 males and 150 females). In this work, we report allele frequencies for the ten X-STR in the Lima population and forensic statistical parameters. Also, the database obtained for the 10 X-STR was compared with data on 15 Iberian and Latin American populations. The genetic diversity analysis showed significant differences in 14 cities evaluated standing it in a single group.