

COMPARISON OF MICROVARIANTS AT DYS458 IN IRAQ, AFGHANISTAN, AND U.S. POPULATIONS

Tabitha Bullock
BAE Systems, 16541 Commerce Dr., King George, VA 22485

DYS458 is one of sixteen commonly examined four basepair Y-STR loci. This poster presents data investigating the occurrence of the microvariants 17.2, 18.2, 19.2, and 20.2 in Iraq, Afghanistan, and U.S. populations. Data from Iraq and Afghanistan was established from profiles generated from detainees and enemies killed in action. Data for the U.S. populations was gathered through published U.S. databases. An elevated occurrence of one or more of these microvariants in the Iraq population is observed when compared to the U.S. and Afghanistan populations.