White-tailed deer DNA links hunter to homicide

Christopher J. Basten

The investigation of a 1997 homicide of a deer hunter near Uniontown, Pennsylvania produced biological evidence from at least three white-tailed deer. A search of a suspect's home produced fresh venison chops and steaks. DNA typing using seven STR loci revealed a genetic match between the steaks and chops and one set of deer remains found at the crime scene. This presentation will compare and contrast the statistical and population genetic issues of human and animal genetic markers for forensic uses.