

Overview of the FBI's CODIS Program

Douglas R. Hares, Ph.D., NDIS Custodian

Federal Bureau of Investigation, Laboratory Division, Quantico, Virginia, USA

The Combined DNA Index System, or CODIS, has matured into a program supporting DNA databases in law enforcement laboratories across the United States. Internationally, there are also 81 CODIS sites in 46 countries. As DNA database programs evolve, CODIS software has and will continue to accommodate the changing needs of the DNA community.

A major component of the CODIS program is operation of the National DNA Index System, or NDIS. NDIS is the National DNA Database. This national index is made up of over 190 DNA laboratories from all 50 States as well as the Department of Justice, the District of Columbia, Puerto Rico, and United States Army Crime Laboratory. Currently there are over 12 million offender DNA profiles and over 500,000 crime scene DNA profiles stored in NDIS.

An update will be provided on the status of the National DNA Index System and the progress of the consortium validation effort for expanding the CODIS core loci in the United States.