

### ***CODIS and NDIS Update***

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 98 CODIS sites in 57 countries. As DNA database programs evolve, the 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 also referred to as the National DNA Database. This national index is made up of over 200 DNA laboratories from all 50 States as well as the Department of Justice, the District of Columbia, Puerto Rico, and the United States Army Crime Laboratory. Currently there are over 16.5 million offender DNA profiles and over 850,000 crime scene DNA profiles stored in NDIS. The use of the CODIS software has resulted in over 400,000 investigations aided.

An update will be provided on the status of the National DNA Index System, future updates to the CODIS software and the FBI's progress toward implementing the Rapid DNA Act of 2017.