STRBASE: 10 YEARS AND BEYOND

John M. Butler

National Institute of Standards and Technology, Biochemical Science Division, Gaithersburg, MD 20899

For more than a decade, the National Institute of Standards and Technology (NIST) has maintained the Short Tandem Repeat DNA Internet Database (STRBase), which is located at <a href="http://www.cstl.nist.gov/biotech/strbase/">http://www.cstl.nist.gov/biotech/strbase/</a>. The purpose of STRBase has been and continues to be an attempt to bring together the abundant literature and information in the human identity testing field in a cohesive fashion to make current and future work easier. New resources that have been recently added will be reviewed.