

EXPERIENCES OF A LARGE DNA DATABANK PROGRAM

Paul B. Ferrara, PhD

Virginia Division of Forensic Science, Richmond, VA



In 1989 and 1990, Virginia passed the first and most comprehensive DNA databank laws in the country. This paper will describe not only the successes of this program but the seemingly endless list of obstacles, issues and decisions that were necessitated as the program evolved. In addition, an analysis of the data which has resulted from this program and the implication of these results on present and future legislation will be discussed.