

CASEWORK EXPERIENCE USING Y-STRs

Cassie Johnson, M.S., Joseph H. Warren, B.A., Judith I. Floyd, B.S., and Rick W. Staub, Ph.D.

Orchid Cellmark, 2600 Stemmons Freeway, Suite 133, Dallas, TX 75207.

Based on several years of experience in offering Y-STR testing at Orchid Cellmark-Dallas, a summary of casework utilizing Y-STRs will be presented to aid analysts, law enforcement personnel, and lawyers in recognizing cases in which Y-STR testing may yield results. This summary will include a description of scenarios in which Y-STR testing is recommended, followed by specific casework examples to illustrate each situation. In addition to discussing the many advantages of Y-STR analysis, the difficulties encountered in Y-STR testing will be addressed. Finally, a strong case will be made for the development of an adjunct Y-STR CODIS database as an investigational aid for law enforcement personnel.