



8th INTERNATIONAL SYMPOSIUM ON HUMAN IDENTIFICATION

September 17-20, 1997

Scottsdale, Arizona

Symposium Topics:

- STRs in Criminal Prosecution
- New Technologies
- Parentage Assessment
- Use of DNA Typing in Forensic Casework
- Population Studies and Databases
- Forensic Applications

Meeting Location:

This is the final year that the Symposium will be held at La Posada Resort in Scottsdale, Arizona. La Posada Resort is located at the base of Camelback Mountain and is a 20-minute drive from Phoenix Sky Harbor International Airport. Conference registration will be accepted by mail, fax or Internet (www.promega.com/genetic/).

Contact: Carol Zabit at (608) 277 2670 (e-mail: czabit@promega.com) or Vicki Schmitz at (608) 277 2639 (e-mail: vschmitz@promega.com)



PRESENTING DNA STATISTICS IN COURT

September 15-16, 1997

Scottsdale, Arizona

A two-day workshop sponsored by Promega Corporation

Instructors: Dr. George Carmody, Carleton University
Dr. Bruce Weir, North Carolina State University

Contact: Carol Zabit at (608) 277 2670 (e-mail: czabit@promega.com) or Vicki Schmitz at (608) 277 2639 (e-mail: vschmitz@promega.com)

Promega Workshops

“DNA TYPING WITH STRS”

FLUORESCENT DETECTION:

August 12-13

November 18-19

December 9-10

SILVER STAIN DETECTION:

August 5-6

November 11-12

Contact: Carol Zabit at (608) 277 2670 (e-mail: czabit@promega.com) for information on Promega's workshops.