

GenePrint™ SHORT TANDEM REPEAT (STR) WORKSHOPS

As part of Promega's worldwide commitment to customer education and exceptional technical support, we are currently offering a series of workshops entitled "DNA Typing With STRs."



Mary Hoenecke demonstrates the proper method of preparing reagents for polyacrylamide gels.

About the Workshops

COMPREHENSIVE

The 2-day workshops provide practical hands-on experience in all aspects of DNA typing using Promega's GenePrint™ STR Systems, including:

- Amplification
- Agarose yield gels
- Polyacrylamide gel preparation
- Polyacrylamide gel electrophoresis
- Silver staining or fluorescent detection
- Data analysis



Workshop participants load their amplified samples onto a polyacrylamide gel for electrophoresis.

WORLDWIDE

GenePrint™ STR workshops are offered throughout the USA and the world.

PROVEN EFFECTIVE

The workshops are well attended and have been extremely popular with our DNA typing customers. Below are comments from former participants.

"The course was extremely well organized. The instructors are excellent and the methodology is elegant...the STR typing method is so well optimized." (Los Angeles Police Department)

"I feel that I can go back to my lab and get reproducible results using these kits...everything worked..." (Biological Laboratory)

"Very well organized!" (Workshop Participant)

"The instructors were very knowledgeable...they were great..." (Texas Department of Public Safety)

HOW TO FIND OUT MORE ABOUT PROMEGA'S WORKSHOPS

- View our Internet site at www.promega.com/genetic/
- Contact Carol Zabit in the USA at
Tel: (608) 277-2670
Fax: (608) 277-2601
e-mail: czabit@promega.com

Promega Symposia and Training Information

LATIN AMERICAN SYMPOSIUM ON HUMAN IDENTIFICATION

August 11-13, 1997

Brazilia, Brazil

STATISTICS WORKSHOP

September 15-16, 1997

Scottsdale, Arizona

8TH INTERNATIONAL SYMPOSIUM ON HUMAN IDENTIFICATION

September 17-20, 1997

Scottsdale, Arizona

2ND EUROPEAN SYMPOSIUM ON HUMAN IDENTIFICATION

June 9-12, 1998

Innsbruck, Austria

Contact: Carol Zabit at (608) 277 2670
(e-mail: czabit@promega.com) for information on any of the symposia listed above.

Promega Workshops

"DNA TYPING WITH STRS"

FLUORESCENT DETECTION:

May 13-14

June 24-25

July 29-30

August 12-13

November 18-19

December 9-10

SILVER STAIN DETECTION:

August 5-6

November 11-12