

Career Opportunity

Promega Corporation, a world leader in supplying biotechnology products for researchers, has an immediate opening for a **European Regional Technical Specialist – Genetic Identity**.

This new position will help Promega achieve sales goals for the Genetic Identity product line within the European region by developing and maintaining relationships with Genetic Identity market leaders and implementing the sales and marketing plan for the region.

The successful applicant will develop and maintain relationships with key contacts, including forensic and paternity laboratories, implement the strategies outlined in the GI sales and marketing plan, work with local branches and distributors to offer technical assistance and represent Promega at various technical sales functions, including tradeshows and symposia.

Qualifications include a broad-based technical background in molecular biology techniques, and DNA typing laboratory experience. A scientific degree in molecular biology, forensic science or equivalent required; advanced scientific degrees preferred. Extensive experience performing DNA typing protocols is required. Proven ability to effectively communicate, in writing and verbally, across all levels of an organization required. English language skills (fluency), both written and spoken required; multiple languages, especially European languages preferred. In order to develop and maintain relationships with key contacts in the industry, regular travel is required. This position will be based in Europe; however, interested candidates should be ready and able to travel up to 50% of the time, primarily throughout Europe.

Promega is a successful growing company, recognized around the world for innovative products and quality customer/technical support. Interested candidates should send a curriculum vitae, with salary requirements, to hr@promega.com or to Promega Corporation, ATTN: Human Resources, 2800 Woods Hollow Road, Madison, WI 53711 USA. For additional information, please visit our website at www.promega.com or www.euro.promega.com.

*Promega Workshops***“DNA TYPING WITH STRS”**

**WINTER/SPRING 1998
MADISON, WI, USA**

FLUORESCENT DETECTION:

February 24-25

March 24-25

April 7-8

SILVER STAIN DETECTION:

March 10-11

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

2nd European Symposium on Human Identification

**Innsbruck, Austria
June 9-12, 1998**

Abstracts for oral presentation
Deadline: March 11, 1998

Contact Vicki Schmitz at (608) 277 2639
(e-mail: vschmitz@promega.com)



Sponsored by
Promega Corporation

