

Unit 5: Solving a Crime Using DNA Analysis

This unit leads students through a crime scene investigation, using RFLP analysis of crime scene sample DNA. Students also have the opportunity to perform basic chemistry analyses as well. For more advanced students who have access to more equipment and reagents, a protocol for STR analysis using polyacrylamide gels is also provided.

Lecture Presentation



Forensic Science DNA Testing

(10 MB)

Laboratory Protocols



Instructor's Lab Manual

(244 KB)



Student's Lab Manual

(110 KB)

Instructor's Guide



Instructor's Guide

(153 KB)