

Tecan's Freedom EVO[®] makes its television debut in *CSI: Crime Scene Investigation*



Sheila Thomas (left) and Barbara Wesley with the Freedom EVO

Freedom EVO for DNA extraction



A Tecan Freedom EVO[®] 150 liquid handling workstation recently starred in a one-hour special episode of the long-running, forensic science-based television show *CSI: Crime Scene Investigation*.

Sheila Thomas, Product Integrator on the hit show that shoots at Universal Studios in California, USA, approached Tecan on behalf of the television program and explained: "It is part of my job to help make the show realistic by finding the correct items for the set. We have a great team of technical advisors who have extensive law enforcement backgrounds but, since technology is constantly changing, we like to stay on the cutting edge by asking for the recommendations of current crime scene investigators who are out in the field every day. We contacted several law enforcement agencies and I also met with various forensic experts at the 2007 International Association of Identifications Conference, who all agreed that Tecan's Freedom EVO platform is one of the best systems available for processing DNA."

Tecan's engineers configured the Freedom EVO workstation for automated DNA extraction from forensic samples and sent the system to Hollywood, where it was installed into *CSI's* on-set laboratory. During the episode, the *CSI* lab technician, Wendy Simms, processes some DNA samples using the Freedom EVO. She has just finished running the samples when special agents Malone, Grissom and Willows enter the lab to discuss the identity of several different cases. The platform is clearly in view throughout the scene and was essential to the storyline because it provided proof that the DNA was not a match for the missing boy.

"I wanted to thank everyone at Tecan for all the hard work they put into getting the Freedom EVO up and running for our show," Sheila said. "Our director is a perfectionist and he was very pleased with how the Freedom EVO performed. We got it in one take!"

The special episode aired in the USA on November 8th 2007.