

## **CALL FOR QUALITY: FORENSIC QUALITY STANDARDS**

Kristina A. Pearson and Kathryn D. Hebble

Promega Corporation, 2800 Woods Hollow Road, Madison, WI 53711, USA

In 2009, the European Network of Forensic Science Institutes (ENFSI), Scientific Working Group on DNA Analysis Methods (SWGDM) and Biology Specialist Advisory Group (BSAG) organizations published a position statement requesting that manufacturers of disposable plasticware and other reagents for the forensic market take precautions to prevent contamination in their manufacturing processes. Here we describe the contamination prevention directives and provide examples of how Promega exceeds all of the guidelines to reduce the chance of human-to-product contamination. Focus is placed on separation of pre- and post-amplification processes, restricted access to labs and QC procedures that ensure lot-to-lot consistency.