

Validation of the Qiagen BioRobot EZ1: An Quick, Automated Method for Extracting DNA.

Kenneth Pfoser, M. Kelly Gannon

Current DNA extraction methods require lengthy extractions and numerous sample manipulations. The Qiagen BioRobot EZ1 is an automated extraction system based on magnetic separation technology requiring minimal "hands-on" preparation utilizing reagent cartridges in an enclosed system. The study evaluated the BioRobot EZ1 for "ease of use", reproducibility, sensitivity, contamination, non-human mixtures, environmental/chemical insults, and non-probative samples.