

A PERFORMANCE EVALUATION AND OPTIMIZATION OF A PROTOTYPE HIGH-THROUGHPUT GDNA EXTRACTION KIT

Caitlin CM Vogelsberg, Crystal M Richt, Jonathan R Galina-Mehlman

Human Origins Genotyping Laboratory, Arizona Research Laboratories, The University of Arizona, Tucson, AZ

High-quality DNA extraction instrumentation and protocols are essential to the consistency and reliability of high-throughput genotyping and analysis services. As a core facility that conducts genotyping on over 100,000 human samples per year, an extraction protocol that produces high quality DNA with minimal handling time is of great importance. We evaluated a novel prototype 96-well vacuum manifold gDNA extraction kit. We compared its DNA isolation performance on 4 different tissue types (bovine bone, liver, blood and human buccal swabs) to that of magnetic particle-based chemistry on a robotic automation platform following the protocols provided by the manufacturers. In addition we optimized the prototype extraction protocol using liver and buccal samples, and evaluated it in parallel with our existing optimized extraction platform. To verify that the vacuum isolation kit produced DNA of suitable quality for PCR amplification, we performed the PCRs on the buccal swab isolations from both methods on one of our laboratory's multiplex PCRs. We visualized the PCR products via fragment analysis on an AB3730, and evaluated for peak quality, peak height and extraneous "noise" or false peaks.

Citations:

Alonso, A., S. Andelinovic, P. Martin, D. Sutlovic, I. Erceg, E. Huffine, L. Fernandez de Simon, D. Albarran, M. Definis-Govanovic, A. Fernandez-Rodriguez, P. Garcial, I. Drmic, B. Rezig, S. Kuret, M. Sancho, D. Primorac 2001. DNA Typing from Skeletal Remains: Evaluation of Multiplex and Megaplex STR Systems on DNA Isolated from Bone and Teeth Samples. *The Croatian Medical Journal*. 42:260-262.

Fahle, G.A. and S.H. Fischer 2000. Comparison of Six Commercial DNA Extraction Kits for Recovery of Cytomegalovirus DNA from Spiked Human Specimens. *Journal of Clinical Microbiology* 38(10): 3860–3863.

Schmerer, W.M., Susanne Hummel, Bernd Herrmann, 1999. Optimized DNA extraction to improve reproducibility of short tandem repeat genotyping with highly degraded DNA as target. *Historical Anthropology and Human Ecology* 8:1712-1716.

Termine, J.D., A.B. Belcourt, K.M. Conn, H.K. Kleinman 1981. Mineral and Collagen-binding Proteins of Fetal Calf Bone. *Journal of Biological Chemistry* 256:10403-10408.

Frame, J and N. Bryant 2007. Optimization of a Qiagen Blood DNA Extraction Kit for use with human buccal swab samples. 18th International Symposium on Human Identification, Poster.