

POWERPLEX® Y 23 PROTOTYPE KIT: AN INTERNAL VALIDATION

Mary Acosta Loyo, Irma Pereira Silva
UEGF, Venezuelan Institute of Scientific Research (IVIC)

The PowerPlex® Y23 prototype kit, was tested with different types of samples, directly or indirectly and using different methods for extracting DNA, such as, salting out (Miller et al.), minikit (QIAGEN) and FTA. For the PCR reaction and the runs on the sequencer, the protocols used were those recommended by Promega with partial modifications in some cases. The following samples were selected for testing: FTA (punch 0.2mm), paraffin embedded biopsies and bone samples. We obtained successful results in at least one of each type of sample. Also, using the same quantity of DNA (0.5 ng), a greater number of systems could be analyzed with the PowerPlex® Y23 with others similar kits. The kit is an excellent tool to the field of human identification. ☺