

GUIDELINES FOR USE OF POWERPLEX KITS ON APPLIED BIOSYSTEMS 3500/3500XL GENETIC ANALYZER

Anupama Gopalakrishnan, Dawn Rabbach, Terri Sundquist, David Yoder and Benjamin Krenke
Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711 USA

Promega's PowerPlex chemistry kits are compatible with many capillary electrophoresis instruments and have been shown to provide reliable and concordant data for human identification.

We provide guidelines for use of the PowerPlex kits on the 3500 instrument for the following:

- generating a matrix for spectral calibration for 4 or 5-dye chemistry
- protocols for running the PowerPlex chemistry kits
- panels, bin and stutter files for data analyses on the GeneMapper IDX software
- global technical support for set-up, validation and troubleshooting of PowerPlex kits