

IMPROVEMENTS TO THE 4-DYE AND 5-DYE MATRIX STANDARDS AND INTERNAL LANE STANDARD FOR 5-DYE POWERPLEX® SYSTEMS

Dawn Rabbach, Mary Baltos, Irina Vyazunova

Promega Corporation, Madison, WI, USA

A new user-friendly version of the PowerPlex® 4C Matrix Standard and the PowerPlex® 5C Matrix Standard that are intended for use on the 3130 and the 3500 Genetic Analysis Systems will be introduced. This user-friendly design combines all matrix fragments in a single tube requiring a single dilution step.

Along with the new format of PowerPlex® 5C Matrix Standard, the existing Promega PowerPlex® five-color STR systems will be retrofitted with a brighter, more stable Internal Lane Standard. The WEN Internal Lane Standard 500 will be replacing the current orange dye, CC5 Internal Lane Standard 500.

Apart from higher signal and stability, WEN Internal Lane Standard 500 has a recommended storage temperature at +4C following initial use. Thereby negating the necessity for thawing the Internal Lane Standard prior to each use.