

Quantifiler® DNA Quantification Kits

Quantifiler® Duo DNA Quantification Kit - P/N 4387746

Developmental validation experiments were performed by Applied Biosystems to evaluate the performance of the Quantifiler® Duo DNA Quantification Kit.

These experiments were performed according to the DNA Advisory Board (DAB) Quality Assurance Standards (October 1, 1998) and the Scientific Working Group on DNA Analysis Methods (SWGDM, July 10, 2003).

Results from these instrument-reagent developmental validation studies have been demonstrated to be fit for purpose in Human Identification testing for databasing, casework, and paternity applications.

Please note the customers should perform their own internal validation studies to establish interpretation criteria and demonstrate that the Quantifiler® Duo DNA Quantification Kit is appropriate and fit for their forensic uses.

