HOW TO IMPROVE STR ANALYSIS USING A NOVEL QUANTIFICATION TECHNOLOGY: MORE THAN DNA QUANTIFICATION

Di Pasquale, F., Cornelius, S., König, M., <u>Bochmann, L.</u>, Prochnow, A., Schnibbe, T., and Engel, H.

QIAGEN GmbH, Hilden, Germany

Quantification in forensic casework analysis is typically the only pre-STR analysis step and could be seen as a quality control step rather than just a quantification of the DNA concentration.

Thus, we present data from our novel human DNA quantification assay product line, called Investigator® Quantiplex Kits, that is available as human (Quantiplex Kit) and human/male (Quantiplex HYres Kit) configuration. Both provide fast and accurate quantification of DNA in forensic database and casework samples with a very high sensitivity and accuracy due to the trusted and validated autosomal multi-copy target 4NS1C and the validated multi-copy target on the Y-chromosome. Detection of inhibitors is ensured by a balanced internal amplification control, without significantly affecting the DNA quantification results. This novel technique of high quality control standards provides a highly accurate assay for DNA quantification in forensics. The assay ensures an improved correlation to the STR results in comparison to other standard real-time PCR methods.

The Investigator Quantiplex product line utilizes a novel PCR fast-cycling technology as well as Scorpion primers that enable rapid results. Using the Rotor-Gene Q system, quantification can be performed in around 50 minutes. ******