

**PARENTAGE TESTING FOR A CHIMPANZEE FAMILY USING ABI AMPFLSTR®-
IDENTIFILER AND ABI AMPFLSTR®- Y-FILER**

Maria Lourdes Andres, Eileen Wu and Patricia Stapleton

DNA Diagnostics Ltd, Auckland, New Zealand

DNA Diagnostics is a human parentage testing lab based in NZ. Recently, Wellington Zoo needed a paternity test done on the newest addition to their chimpanzee family.

Since chimpanzees are the closest living evolutionary relatives to humans there is enough commonality in our DNA to warrant the use of ABI AmpFISTR® Identifiler (Identifiler) and ABI AmpFISTR® Y-Filer (Y-Filer) to determine the paternity of the baby.

DNA was extracted from the hair samples provided from the three possible fathers, mother and male baby. In Identifiler, amplification products from eight regions of DNA showed variation between the five chimpanzees. In Y-Filer, the males showed one amplification product at each of the seven regions and unexpectedly the female showed amplification at two regions.

Based on the Identifiler data, one of the potential fathers could not be excluded from paternity, but a complication arose in the Y-Filer data where he did not match the baby at one locus. We believe this to be a mutation.