

ANALYSIS OF AUTOSOMAL AND Y-CHROMOSOME STR FROM CIGARETTES BUTTS AFTER TREATMENT WITH FINGERPRINT REAGENTS

Reena Roy, PhD

St. Louis County Police Crime Lab, Clayton, MO



Cigarettes smoked by a man and a woman were stored in paper envelopes for a week and then fingerprinted using ninhydrin and Super Glue[®] fuming. DNA was successfully extracted from approximately 5mm paper of the butt end of each cigarette using the DNA IQ[™] System. Purified DNA was amplified using the PowerPlex[®] 16 System for the detection of the autosomal STR. DNA extracted from the cigarette butts smoked by the male individual was also subjected to amplification by the Y-PLEX[™] 6 and Y-PLEX[™] 5 amplification kits.

Complete autosomal and Y-STR profiles were obtained from the cigarette butts after treatment with the fingerprint reagents.