

FLUORESCENT FOUR-STATE MSI MVR-PCR MAPPING OF DNA FROM CHEWING-GUM IN CASEWORK

Peter Hau¹ and Nigel Watson²

¹*Lothian and Borders Police Forensic Science Laboratory, Edinburgh, Scotland*

²*Forensic Science Unit, Strathclyde University, Glasgow, Scotland*



Minisatellite Variant Repeat Mapping by PCR for the MS32 and MS1 loci have been described previously (Jeffreys *et al*, 1991 and Hau & Watson, 2000). The four state mapping method makes use of four repeat specific PCR primers directed to different repeat sequence types. Therefore a high degree of variation in the minisatellite is revealed. MS1 is the most polymorphic locus used in forensic science. This poster reports the application to the DNA recovered from discarded chewing-gum found at the scene of a crime.