

TOUCH DNA - THE FORENSIC USE OF SOLID AND POROUS MATRICES COATED WITH POLYANILINES IN THE SELECTIVE SEPARATION OF BIOPOLYMERS FROM NUCLEIC ACIDS.

Belen Rodriguez Cardozo¹, Stella Maris Jaureguiberry² and Gabriel Oscar Boselli³

¹Banco Nacional de Datos Genéticos, Hospital Dr. C.G. Durand, J.B. Ambrosetti 743 – PB, 1405 Ciudad de Buenos Aires, Argentina

²Superintendencia de Policía Científica, Laboratorio de Genética Forense, Ministerio de Seguridad Policía de la Provincia de Buenos Aires, Calle 91 entre 48 y 46 nro 1829 2do piso, 1653 San Martín, Provincia de Buenos Aires, Argentina

³BioSystems S.A, Dorrego 673, 1414 Ciudad de Buenos Aires

Selecting appropriate methods of collection and extraction of trace DNA samples from a crime scene could be relevant during the investigative process.

Actual extraction methods are based on DNA binding using different sorbents, which could cause significant loss of the starting material.

In the present work we performed comparative studies of a new non-binding separation method (Fingerprint DNA Finder Kit – FDF® Kit) that uses porous matrices associated with polyanilines nano-layers, which are able to retain selectively biopolymers and potential PCR inhibiting substances, while nucleic acids are never bound and remain in solution, collecting all DNA from nucleated cells, keratinocytes and even cell-free nucleic acids (basically free-floating DNA fragments not encapsulated in the cell nucleus).

Dealing with comparative studies of trace lift protocols implies the use of controlled samples. The preparation of control samples using dilutions of blood or saliva, and applied in liquid drops over different surfaces, does not reproduce exactly the conditions of epithelial cells or free DNA contained in the sweat. We use here real touched or manipulated objects as the only DNA source (touch DNA).

The FDF® Kit demonstrated being more effective, not only in DNA yield, but also in the quality of the profiles, which were always clear and devoid of artifacts. The use of this new kit associated with adhesive trace lifters was sensitive enough to be applied in real casework. **⌘**