

INTERPOL's DNA Unit: Tools and Services for International DNA Data Exchange

Andrea Fischer, PhD and Susan Hitchin, PhD

INTERPOL (International Criminal Police Organisation)

INTERPOL, as the world's largest international law enforcement organisation, assists police in its 190 member countries in the fight against international crime by making available tools and services for the international exchange of police data including forensic DNA data.

Tools and services for DNA data exchange range from providing a central DNA database, hosting bilateral DNA data exchanges, to contributing to the pursuit of wildlife poachers through DNA-based methods on animal remains.

INTERPOL's DNA database, maintained by the DNA unit in Lyon, France, is used to link transnational perpetrators to crimes committed in other countries as well as providing a connection between DNA profiles of unidentified human remains and missing persons on an international scale. Additionally, INTERPOL is expanding its DNA capabilities to include kinship matching that can provide further support for international missing persons investigations.