

SAO PAULO STATE UNIVERSITY AND PUBLIC DEFENDER: TWENTY NINE MONTHS OF PATERNITY CASES HELPING CHILDREN OR ADULTS WITHOUT FINANCIAL SUPPORT FOR PAYING THE TEST

Regina Maria Barreto Cicarelli, Fernanda Silva Polverari, Isabela Brunelli Ambrósio, Simone Busko Valim

São Paulo State University – UNESP Laboratório de Investigação de Paternidade - Núcleo de Atendimento a Comunidade - Faculdade de Ciências Farmacêuticas - Universidade Estadual Paulista

The UNESP Paternity Research Laboratory has an agreement with the Public Defender's Office of the State of São Paulo and since 2015, attending to the population that is unable to solve cases of affiliation: trios, duos, reconstructions and exhumations. The tests have been carried out with Promega kits: PowerPlex21, PowerPlex Fusion and, more recently, PowerPlex Fusion 6C. Exclusion confirmations are analyzed by sex chromosomes (PowerPlex Y23 and Decaplex X-STRs). During this period, 1283 cases (including few private ones) and 24 post-mortem cases were collected, with 1383 reports being prepared and 4608 individuals analyzed. The percentages of inclusion were always higher than those of exclusion and/or inconclusive. Data concerning the question of whether the mother informs or not if there is another biological father will also be discussed.

Financial support (partial): FAPESP (2016/10138-8)