
CONQUERING MT. EVEREST: COMPLETING THE ALL-FELON OFFENDER BACKLOG

Robin L. Rothove

Missouri State Highway Patrol Crime Laboratory, 1510 East Elm Street, Jefferson City, MO 65101

The All Felon law was implemented in Missouri in January 2005 and created an immediate backlog of a 100,000 samples. The program rapidly evolved from analyzing 2,500 samples each year to 4,500 samples each month. Discover how the Missouri State Highway Patrol completed the backlog within two years without robotics or outsourcing. Learn how the MSHP completed the backlog within two years without robotics or outsourcing! Discover how the procedure, budget, equipment and employees made it all possible while also uncovering the “dos” and “don’ts” associated with completing a seven-year project in two years.