
The Innocence Project- An Update

Maddy deLone
Executive Director of the Innocence Project

The Innocence Project is a nonprofit legal clinic that uses DNA evidence to exonerate wrongfully convicted individuals, will discuss the legal and policy issues related to post-conviction DNA access. After delivering a brief overview of the Project, she will discuss the 207 exonerations (to date) that relied upon DNA evidence. She will outline the causes of these wrongful convictions and then move into a policy discussion of DNA access laws in the United States, with a focus on the states where there are weak laws or no laws at all. Her presentation will conclude with a question-and-answer session.