The Killer of Little Shepherds

Douglas Starr

Boston University, Boston, Massachusetts, USA

At the end of the nineteenth century, serial murderer Joseph Vacher, known and feared as "The Killer of Little Shepherds," terrorized the French countryside. He eluded authorities for years—until he ran up against prosecutor Emile Fourquet and Dr. Alexandre Lacassagne, the era's most renowned criminologist. The two men, intelligent and bold, typified the Belle Époque, a period of immense scientific achievement and fascination with science's promise to reveal the secrets of the human condition.

In this presentation, Douglas Starr will revisit Vacher's infamous crime wave, interweaving the story of how Lacassagne and his colleagues were developing forensic science as we know it. We see one of the earliest uses of criminal profiling, as Fourquet painstakingly collects eyewitness accounts and constructs a map of Vacher's crimes.