

A COLD CASE REVIEW INTO THE MIND OF SERIAL KILLER JOSEPH NASO

Dawn Romano, Contra Costa County Crime Lab

In 2010 a search of Joseph Naso's house as a result of a probation violation led to the gruesome discovery of a treasure trove of photographs of young women appearing drugged or dead, a "rape" diary, and a top ten list. An investigation conducted on Naso's top ten list uncovered a link to the murder of a series of young women in northern California over a span of several decades. This talk will outline the entire investigation from probation violation to Naso's eventual conviction for the rape/murder of four of his victims.