

MUHAMMAD AND MALVO: THE SERIAL SNIPER CASE OF OCTOBER 2002

Brendan Shea, M.S., and Richard Guerrieri, M.S.

Federal Bureau of Investigation Laboratory Division, Quantico, Virginia 22135, USA

The serial sniper attacks that occurred in October 2002 in Maryland, Virginia, and the District of Columbia caused region-wide levels of fear and tension and effected the lives of millions of residents in the area. This presentation will provide insight into the sequence of events surrounding the sniper attacks, some of the contacts that took place between the perpetrators of these acts and law enforcement, and discuss some of the results obtained by the FBI Laboratory. Specific emphasis will be placed on the DNA examinations conducted and the unique items and examinations that yielded probative DNA typing results.