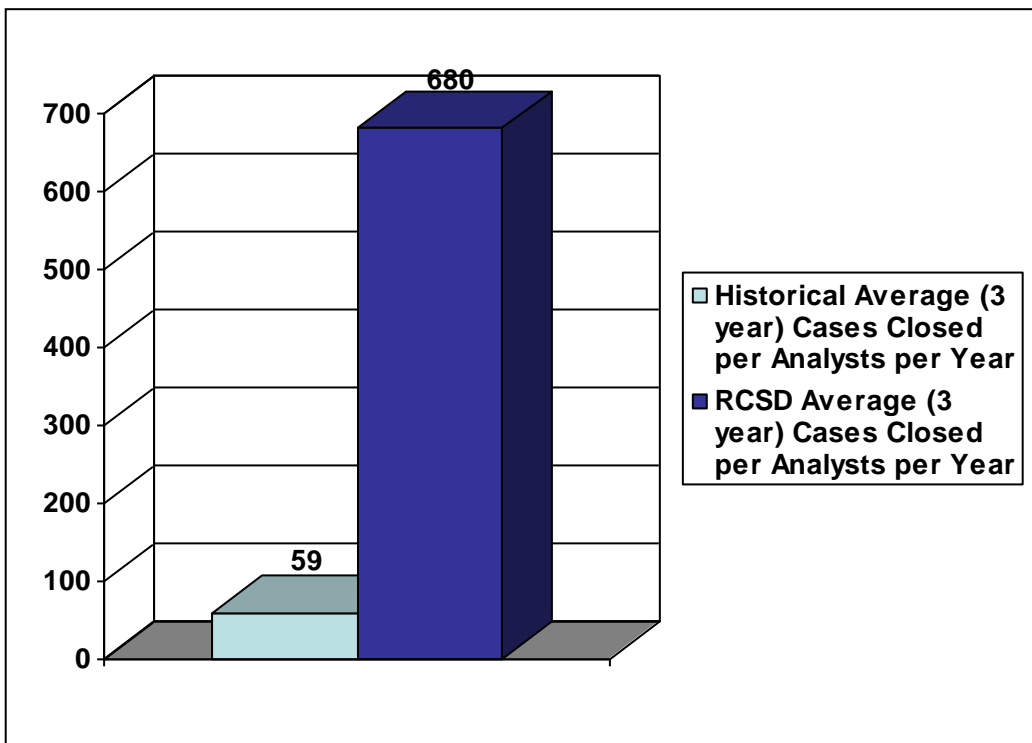


## PROCESS ENHANCEMENTS AT RICHLAND COUNTRY SHERIFF'S DEPARTMENT

Grayson Amick  
Richland County Sheriff's Department

The Richland County Sheriff's Department's DNA laboratory has implemented process enhancements over its 5 year existence that has resulted in an improvement in casework efficiency. A three year average of cases completed exhibited an 11.5-fold increase versus historical data. Hence an average of approximately 680 cases was completed per analysts per year compared to an average of approximately 59 cases per analyst per year. In addition, analysis time and cost have decrease approximately 8-fold and 2-fold respectively over a 5 year period. These improvements are a result of process improvements, integration of automation, technology advancements, and novel reporting conventions.



2009, RCSD, Gray Amick