LEVERAGING LABORATORY DATA IN DECISION MAKING

Robin Gall, and Cynthia Shannon

Abbott Informatics

Laborato	pries have mountains and mountains of data, but is it being effectively leveraged i
decision making? This is one challenge in the forensic lab today. Have you asked:	
	How do I gather relevant insights from this data?
	Am I leveraging all of the available data in this decision?
	What is causing the spike in turnaround times?
	Where is the bottleneck in the process?
	What resources are needed, personnel, equipment, reagents, all of them?

You have decisions to make. Do you have the necessary information to make the best decision? How many systems do you currently use to extrapolate data to make an informed decision?

Advanced Analytics is an integrated solution with STARLIMS that provides graphical and numerical representations of your data. AA allows you to dynamically interact with the information. For example, do you want to show how grant funds were utilized in your laboratory? Advanced Analytics will display the resources associated with a grant. Want to know what cases were analyzed using those resources? Simply click on the resource and the cases and exhibits associated with the resource are displayed. Furthermore, you may export the information in a graphical display, pdf, or Excel, so it may be inserted into a report.

Advanced Analytics provides new insights, so you can draw meaningful conclusions and make better decisions. You can analyze laboratory operations in new ways; identify problem areas, drill down to find underlying root cause and start to formulate solutions; identify trends and patterns over time; and increase productivity and laboratory efficiency.

Advanced Analytics and STARLIMS allows you to answer tough questions with real time information; conduct root cause analysis and quickly understand the impact of an instrument, reagent, or analyst issues. With Advanced Analytics the opportunity to mine your data is endless, because you have the ability to build new workbooks to understand your data. With the ability to have meaningful information, you can focus on the important stuff!