



Tamine Hosseini (third from left) and colleagues in the Department of Clinical Chemistry at the Sahlgrenska University Hospital

Automated pre-analytical blood sample sorting

The Sahlgrenska University Hospital in Gothenburg – one of the largest hospitals in Sweden – is saving considerable time using a Genesis FE500™ workstation for the pre-analytical processing of blood samples.

At Sahlgrenska, the sample reception section of the Department of Clinical Chemistry is responsible for sorting and processing all hospital samples that arrive for laboratory analyses, including microbiology, immunology and serology, as well as the clinical chemistry samples. The laboratory's Genesis FE500 workstation processes EDTA and citrate coagulation tubes for the full range of departments, as well as preparing some samples for archiving, automating the centrifugation and decapping of the sample tubes in preparation for subsequent analysis.

“Before we had the Tecan system, we had to process these samples manually,” explained Tamine Hosseini, head of sample reception. “I would estimate that we are now saving around 30 to 40 per cent of the time originally spent in sorting, centrifugation and decapping.

The sample barcodes are scanned and registered into our laboratory information system (LIS) along with all the patient information and analysis requests. Once loaded onto the workstation, the barcode specifies the tests needed for each tube and the samples are sorted and processed accordingly.”

“Our staff enjoys working with the Genesis FE500 workstation, largely because it needs virtually no daily, and only minimal weekly, maintenance work, compared to the analytical instruments, which require frequent calibration and extensive maintenance routines,” Tamine added.

“We have had the Genesis FE500 system for one year and plans are now underway to use it for processing urine samples in the near future to make the most of the capabilities this instrument offers.”

For more information on Tecan's Genesis FE500, visit www.tecan.com/genesisfe500