



# Benefits of GSP + Specimen Gate Lab

GSP Workstation is an instrument-level software product intended to assist with the usual tasks associated with using the GSP in a laboratory. Revvity's Specimen Gate® (SG) software is designed to simplify the data collection and workflow processes involved with receiving specimens, screening for abnormalities, managing patient information, generating specimen reports, and following up abnormal and unsatisfactory specimens. Some of the benefits of pairing the GSP instrument with Specimen Gate® Laboratory software are listed below:

Feature	GSP WS instrument software	GSP with SG lab
Instrument connections	Limited to 1 GSP	Many; includes non-GSP as well
Smart positive ID™	NO	YES
Fixed cutoffs	YES	YES
Percentile cutoffs	NO	YES
Demographics based cutoffs (Requires Evoya)	NO	YES
Specimen details in history window	NO	YES
Open another assay from history window	NO	YES
Supported acceptance steps	Calculated, Accepted	Calculated, Inspect, Checked, Accepted
Cutoff analyzer	NO	YES
Initial samples in RV request list	NO	YES



- Smart Positive ID enables the software and punchers to communicate. Mistakes can be corrected during the punching process, when the GSP is interfaced with Specimen Gate® Lab.
- GSP Workstation is intended for use with a single instrument. As a Total Process Management Solution, Specimen Gate® Laboratory is intended to address a wide range of instrumentation and connectivity needs.
- Specimen Gate® Laboratory provides a wide range of configuration options to allow customer-specific configuration.
- Revvity now offers a cloud based solution with Evoya that can provide additional downstream supporting features.