



Clinical Diagnostics Overview

Summary	Bio-Rad's Clinical Diagnostics Group develops, manufactures, sells, and supports a large portfolio of products for medical screening and diagnostics. Bio-Rad is a leading specialty diagnostic company, and its products are recognized as the gold standard for diabetes monitoring and quality control (QC) systems. The company is also well known for its blood virus testing and detection, blood typing, autoimmune and genetic disorders testing, and internet-based software products.
Customers	Bio-Rad's primary clinical diagnostics customer base consists of hospital laboratories, reference laboratories, physician's laboratories, government agencies, and other diagnostics manufacturers.
Products	<p>The company's clinical diagnostics products incorporate a broad range of technologies used to detect, identify, and quantify substances in bodily fluids and tissues. The results are used as aids to support medical diagnosis, detection, evaluation, and the monitoring and treatment of diseases and other medical conditions.</p> <p>Quality Control Products Bio-Rad offers the world's most comprehensive menu of quality control products, the leading QC data management software, and unparalleled customer services.</p> <p>Autoimmune Testing Bio-Rad is a leading supplier of diagnostic products for autoimmune testing, helping clinical laboratories and doctors provide patients with more accurate diagnoses and faster, more effective treatment solutions.</p> <p>Blood Virus Testing Bio-Rad is one of the top three suppliers of testing systems for HIV and hepatitis for use by hospitals and blood banks throughout the world.</p> <p>Diabetes Monitoring Bio-Rad is the number one supplier of the glycosylated hemoglobin test that is recommended quarterly for those diagnosed with diabetes. Bio-Rad's sophisticated systems and technical support enable doctors to monitor a patient's status, and ultimately to help patients in their fight to effectively manage and control this disease.</p> <p>Genetic Disorders Screening Bio-Rad plays an important role in screening newborns as well as expectant parents for genetic diseases. Bio-Rad's genetic disorders product line includes a highly specialized line of assays, software, and automation.</p> <p>Immunohematology Screening To improve blood typing reliability and efficiency in private laboratories, hospitals, and blood banks world wide, Bio-Rad offers the gold standard in immunohematology reagents and instrumentation.</p> <p>Microbiology Testing Bio-Rad draws on more than 100 years of experience and expertise through its partnership with the Institute Pasteur in Paris. It offers a full range of sophisticated, reliable, and cost-effective tools for the diagnosis of microbial infections—from traditional culture media to innovative rapid diagnostic assays.</p> <p>BioPlex® 2200 System Launched in 2006, the BioPlex 2200 system is the industry's first and only fully automated, random access multiplex testing platform to generate multiple results from a single patient sample. The system offers significant laboratory cost savings through platform consolidation, decreased labor, and reduced reagent consumption.</p>