



The Universal Electrophoresis Gauge can be used to measure the dimensions of various components used in electrophoresis equipment, such as spacers, combs, and glass tubes.

To use, insert the part to be measured into the appropriate slot on the gauge and move it to the left until it stops. This will give you the approximate dimension of the part. The part should be held perpendicular to the gauge and not forced so that the gauge bends.

The Universal Electrophoresis Gauge also has a centimeter scale/ruler to be used for general measuring purposes.

Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, CA 94547