

## Model RGC-100 and RGC-150 Series Digital Vacuum Gauge Quick Start Guide

The Varian RGC-100 and RGC-150 series gauges are compact digital vacuum sensing instruments that use a thermocouple gauge tube to sense vacuum and display the reading in available units.

1. Verify you have received all components:
  - Controller with LED display
  - Cable that connects display controller and 531 Sensor
  - Varian type 531 Sensor
  - Universal Plug - Power supply
2. Plug your display controller into wall power.
3. Verify display lights up.
4. Plumb your VARIAN 531 gauge tube into the vacuum manifold with the pipe threads facing down, and the octal pins facing up.
5. Connect your gauge tube cable to the display controller on one end, and to the gauge tube with the octal connector on the other end
6. Verify appropriate readings. For example, a chamber exposed to atmosphere might have a reading of 600-760 Torr, which is appropriate.

For the full instruction manual, please see the latest documentation at: [www.varianinc.com/vacuum](http://www.varianinc.com/vacuum). Choose *Downloads*, then *Vacuum Measurement Instrumentation*.

