

# DVG50

## Digital Vacuum Gauge

# INSTRUCTION MANUAL

---

### INSTALLATION

Please read the following instructions thoroughly before attempting to install the instrument.

### RECEIVING

Inspect the Digital Vacuum Gauge for any obvious signs of damage due to shipment while unpacking. Immediately advise the delivering carrier if any damage is suspected.

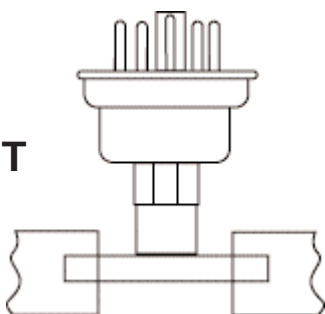
### POWER REQUIREMENTS

The Digital Vacuum Gauge operates on 15 VDC, 400 mA. Models shipped within North America are accompanied with a 115VAC, 60 Hz power transformer. In all other areas, the appropriate transformer suited to the area power requirements is required.

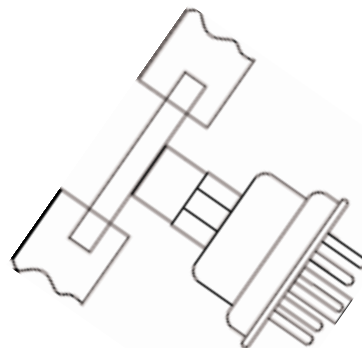
### GAUGE TUBE INSTALLATION

When mounting the gauge tube in the vacuum system, the preferred orientation is a vertical position, to prevent accumulation of liquid and solid particles.

**CORRECT**



**INCORRECT**



The gauge tube is installed in the vacuum system by cutting the vacuum hose and inserting the "T" connection. The connection should be made vacuum-tight by fastening with hose clamps.

### OPERATION

#### READING THE VACUUM GAUGE

When the vacuum level is between 50 and 1 Torr and a decimal point is showing, vacuum is read in units of Torr. If no decimal is seen, vacuum is read in mTorr or microns.

#### EFFECT OF THERMAL CONDUCTIVITY

The Digital Vacuum Gauge is originally calibrated in dry air. This calibration is a function of thermal conductivity. Solvent vapors have different thermal conductivities and will provide readings with a slight variance when compared to absolute pressure. For most application involving the Digital Vacuum Gauge, the vacuum levels noted on the gauge will serve as a relative measure and not absolute.