

# 9964-033

## CO<sub>2</sub> Tank Connector Block

### Installation Instructions for the LI-6400 Portable Photosynthesis System

#### General Description

The 9964-033 CO<sub>2</sub> Tank Connector Block is used with an external CO<sub>2</sub> source and the 6400-01 CO<sub>2</sub> Controller (located in the LI-6400 console) to provide automatic control of CO<sub>2</sub> in the LI-6400. The tank connector block replaces the external CO<sub>2</sub> source assembly, and is useful in situations where a CO<sub>2</sub> source of longer duration than can be provided by the CO<sub>2</sub> mini-cartridges is needed.

The tank connector block is designed for use at pressures between 180 and 220 PSIG of CO<sub>2</sub>. **Do not exceed 250 PSIG CO<sub>2</sub>, as the pressure relief valve may vent.**

The tank mounting block is fitted with a 1/8" male NPT to 1/8" tubing fitting. This fitting is installed with a 10 SCCM flow restrictor. **Do not remove this fitting.** A 1/8" to 4mm compression union is also provided for users who may be unable to obtain 1/8" copper tubing. Directions for installing the tank connector block to a CO<sub>2</sub> source using 4mm copper tubing are given on the reverse.

#### Installation

1. Remove the red protective plastic cap covering the CO<sub>2</sub> inlet on the LI-6400 case, between the CO<sub>2</sub> and H<sub>2</sub>O scrub tubes.
2. The CO<sub>2</sub> tank connector block is mounted on the left side of the LI-6400 console, between the CO<sub>2</sub> and H<sub>2</sub>O scrub tubes. Make sure that the O-ring seal on the back of the block is properly seated. Tighten the two knurled knobs on the mounting block to secure the assembly to the console.
3. Insert a length of 1/8" copper tubing through the 1/8" connector nut and the ferrule (Figure 1). **Important:** Note the orientation of the ferrule. One of the tapered ends of the ferrules is longer than the other; the long end must be oriented toward the connector on the mounting block. When the nut is tightened onto the connector, the ferrule will be permanently crimped to the copper tubing, and you will not be able to remove it.

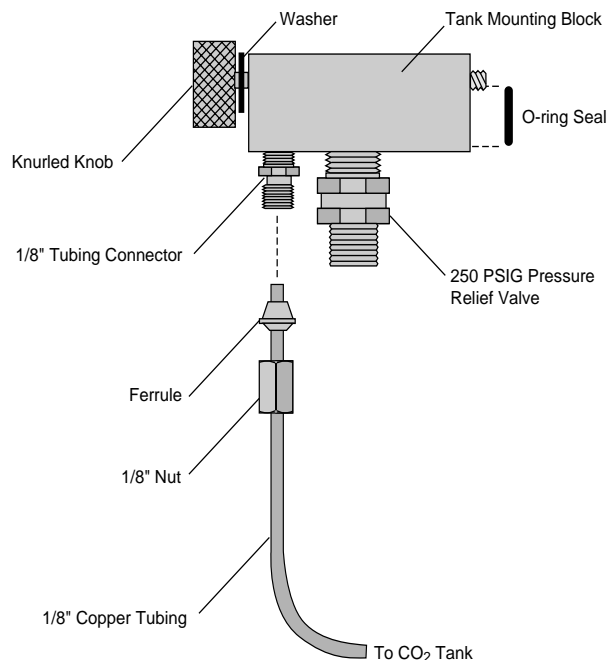


Figure 1. Insert tubing through nut and ferrule. Note proper orientation of ferrule.

4. Tighten the nut until snug, plus 3/4 of a turn.
5. Connect the other end of the copper tubing to your CO<sub>2</sub> source. Adjust the tank pressure to between 180 and 220 PSIG.

## Installation Using 4mm Copper Tubing

If you are unable to obtain 1/8" copper tubing, it is possible to connect the tank connector block to a CO<sub>2</sub> source using 4mm tubing and the compression fitting (LI-COR part #300-04439) included with the tank connector block.

1. Install the tank connector block and the short length of 1/8" copper tubing (included) as described in steps 1-4 above.
2. Use the 1/8" to 4mm compression fitting to connect the 1/8" and 4mm tubing, as shown in Figure 2. Be sure to orient the ferrules correctly; the narrow tapered end of each ferrule must be oriented *toward* the compression fitting. Tighten the nuts on the compression fitting until snug, plus 1 1/4 turn.
3. Connect the 4mm tubing to your CO<sub>2</sub> source. Adjust the regulator pressure to between 180 and 220 PSIG.

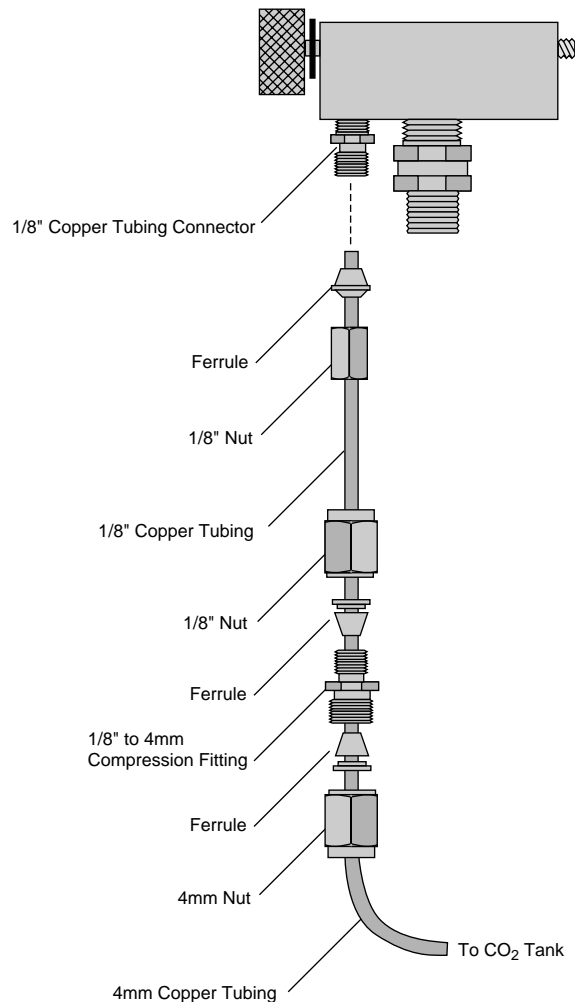


Figure 2. Use the compression fitting to connect 1/8" and 4mm tubing.

**LI-COR®**

LI-COR, inc. ● Environmental Division ● 4421 Superior Street ● P.O. Box 4425 ● Lincoln, Nebraska 68504 USA  
 Phone: 402-467-3576 ● FAX: 402-467-2819  
 Toll-free 1-800-447-3576 (U.S. & Canada)

PPS205