

Setting up your Agilent 6850 GC

1 Remove protective caps from TCD/FID and vents.

2 Connect all cables.

3 Remove plugs and install column (label to front).

4 If available, install cryo cooling coil. Bend bracket to secure coil.

5 If available, install your sampler and cable.

6 Remove caps and connect your gases. If valve installed, connect sample lines.

7 Set gas source pressures and check for leaks.

Gas	Recommended	Maximum
Helium (He)	345 kPa (50 psi)	690 kPa (100 psi)
Hydrogen (H ₂)	345 kPa (50 psi)	690 kPa (100 psi)
Air	345 kPa (50 psi)	690 kPa (100 psi)
Makeup	345 kPa (50 psi)	690 kPa (100 psi)
Actuator air	310 kPa (45 psi)	345 kPa (50 psi)

8 Turn on the GC. If available, connect cryo filter and coolant.

9 Load CHECKOUT method and wait until "Ready..."

10 For manual injection, prepare the checkout sample and start.

11 If available, prepare checkout sample and start.

12 Verify chromatographic peaks.

If available, plug in the control module.

>44 cm (17.3 in)
>21 cm (8 in)
>29.5 cm (12 in)
H: 51 cm (20 in)
W: 29 cm (11.4 in)
34 cm (13.4 in) CO₂
D: 57 cm (22.4 in) G2680
Wt: <23 kg (51 lb)

Instrument Safety Symbols

For more details, consult the information provided on the 6850 User CD-ROM and other accompanying documentation.

www.agilent.com