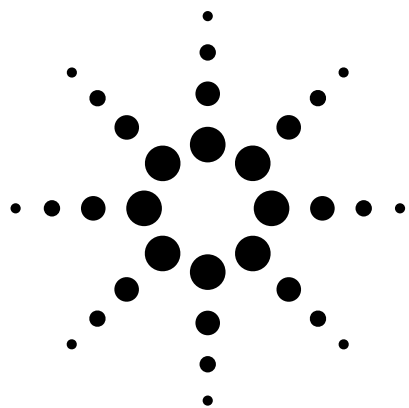




Installing an Auxiliary Pressure Control Module

Agilent 6850 GCs

Accessory G3349B

This kit contains:

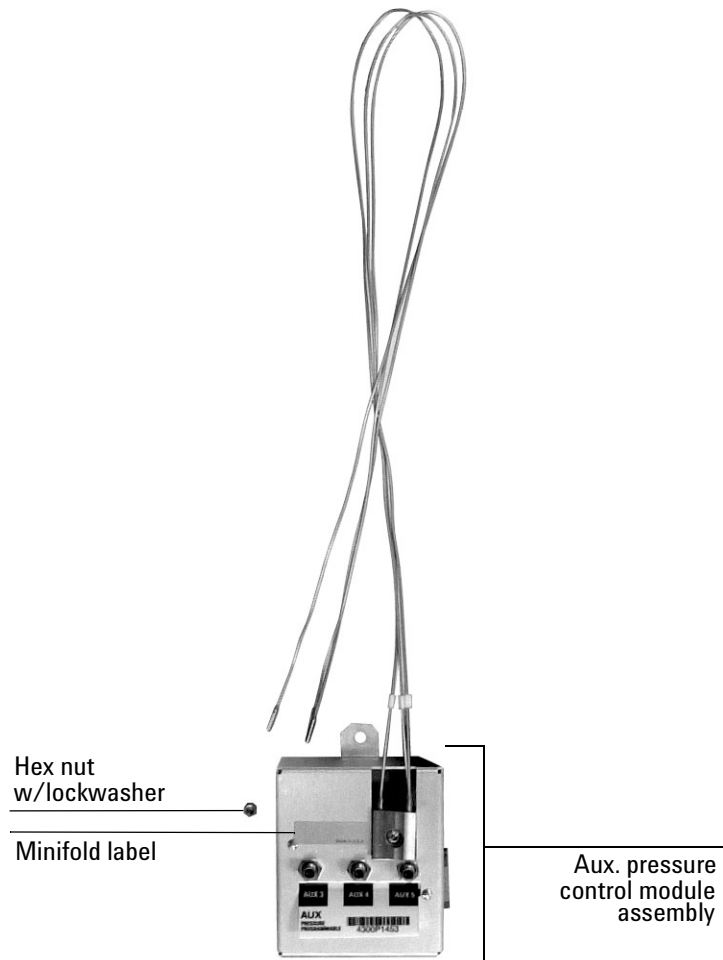
Description	Quantity
Cable tie .062-.625 diameter	2
Hex nut w/lockwasher	1
Disposable wrist strap	1
Minifold label	1
Restrictor kit	1
• FID air restrictor (brown)	3
• Blue dot restrictor	3
• O-rings (6 pk)	1
• Tube zero restrictor	3
• External H ₂ flow limiter	2
• External Flow Restrictor Install Guide	1
Aux. press. control module assembly (with covers)	1
Back panel replacement for the GC	1
6850 GC user information CD	1
ChemStation software CD	1
Installation sheet (this document)	1

The module is factory-assembled. Do not disassemble it during installation.



Agilent Technologies

Parts identification



Required tools

- Torx® T-20 screwdriver
- Needle-nose pliers

Overview

Caution

Before starting, review the safety information listed at the end of this document.

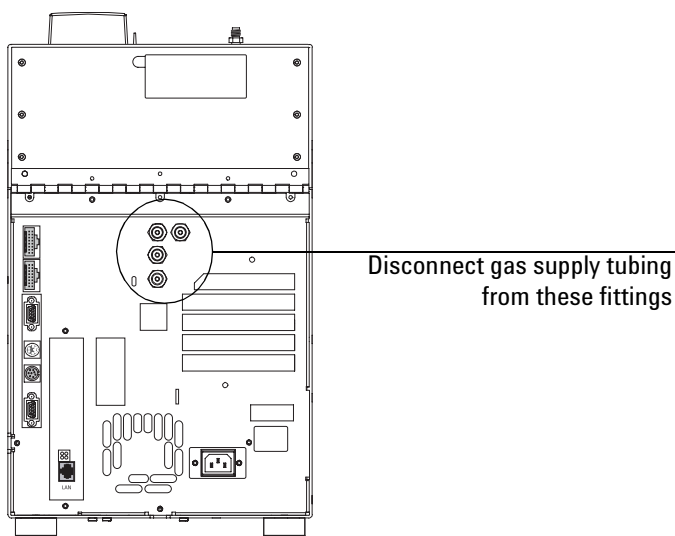
1. Disconnect the GC.
2. Install the new module.
3. Restore the GC to operating condition.

Disconnect the GC

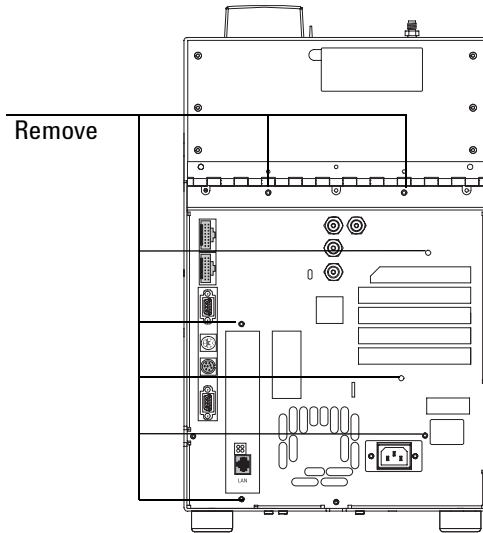
WARNING

Hydrogen gas is flammable and potentially explosive. Before replacing the module, turn off the hydrogen gas at the source.

1. Turn off the GC and unplug the power cord. Allow time for all heated zones to cool.
2. Disconnect all cables from the GC.
3. Turn all gases off at their sources. Disconnect the carrier and detector gas tubing from the back panel of the instrument.



4. Remove the seven T-20 screws on the back panel. Save the screws.



5. Tilt the top of the panel away from the GC.
If you have a Series II GC, save this panel. Otherwise, set this panel aside because you will no longer use it. A new panel comes with the kit.

Install the new module

Caution

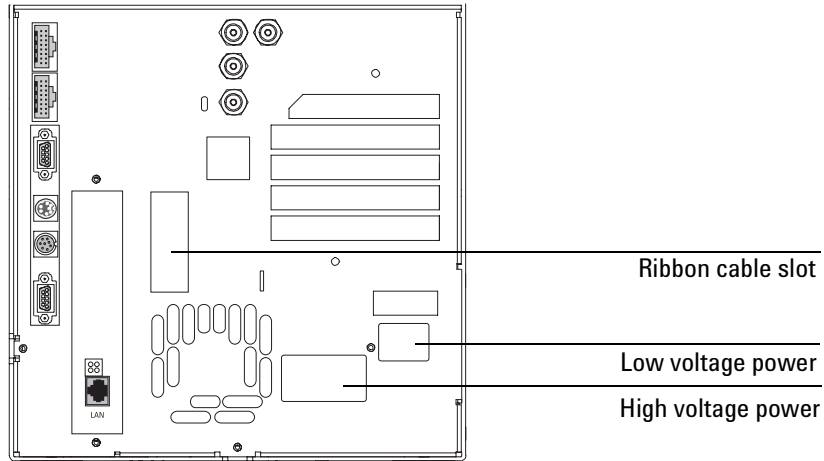
Always hold the module by its support bracket to avoid damaging board components.

1. Change one or more auxiliary channel frits, if needed. See “Changing the flow restrictors (optional)” on page 8.

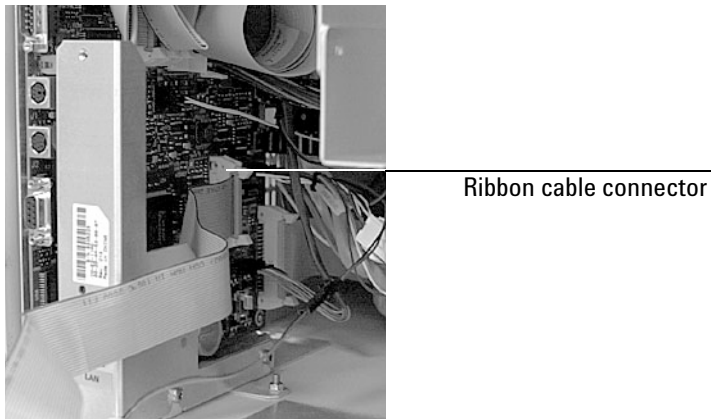
Installing an Auxiliary Pressure Control Module

Install the new module

2. If you will be using the back panel from the kit, use needle-nose pliers to remove the knockouts for the ribbon cable and power cable slots.



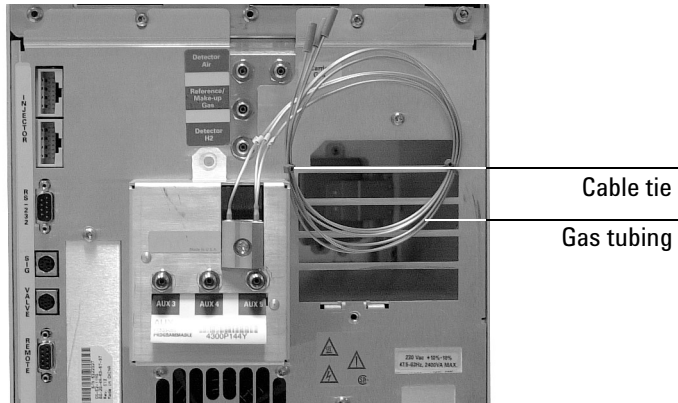
3. Remove the tape from around the new control module and thread the ribbon cable through the slot in the back panel.
4. Attach the ribbon cable to the pneumatics control board (at P9).



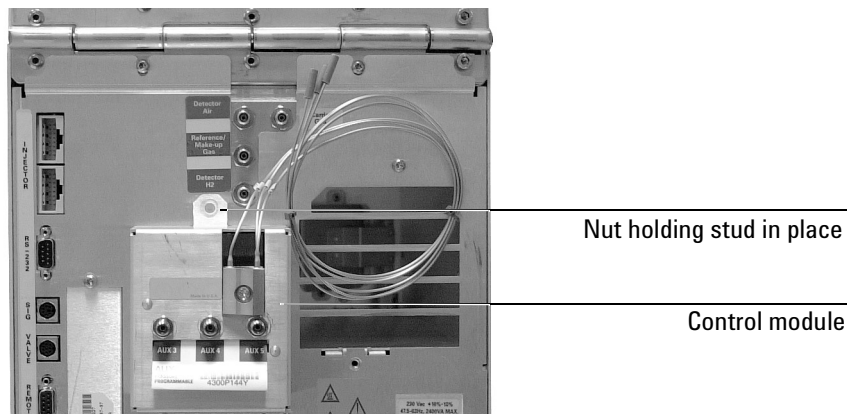
Installing an Auxiliary Pressure Control Module

Install the new module

5. Gather the excess gas tubing from the gang fitting block in a loop and tie it together with a cable tie.



6. Attach the control module to the back panel by hooking the tabs on the module just inside the ribbon cable slot. Position it so you can insert the threaded stud through the hole in the panel. On the inside, use the hex nut with lockwasher to hold the stud in place.
7. Install the back panel.



Changing the flow restrictors (optional)

The auxiliary channels are controlled by a pressure setpoint. To work properly, there must be adequate flow resistance downstream of the pressure sensor. The auxiliary channel pneumatics manifold provides a frit-type restrictor for each channel.

Four frits are available:

Frit marking	Flow resistance	Agilent part number
Blue Dot	High	19234-60660
Red Dot	Medium	19231-60770
Brown Dot	Low	19231-60610
None (brass tube)	Zero	G1570-20540

The red dot frit is in all three channels when the instrument is shipped.

Figure 1 and Figure 2 show approximate pressure/flow relationships for the three dot frits, assuming there is no significant resistance downstream of the frits.

If the Zero resistance frit is installed, the user must provide flow resistance downstream and generate the pressure/flow relationships.

WARNING

When hydrogen is used, dangerously high flows are possible if insufficient flow resistance is provided downstream of the supply tube. Always use either the High (blue dot) or Medium (red dot) frit with hydrogen. You may also want to use the restrictor kit that comes with this accessory. Its installation document describes using it for the 6890 GC, but still applies to the 6850 GC.

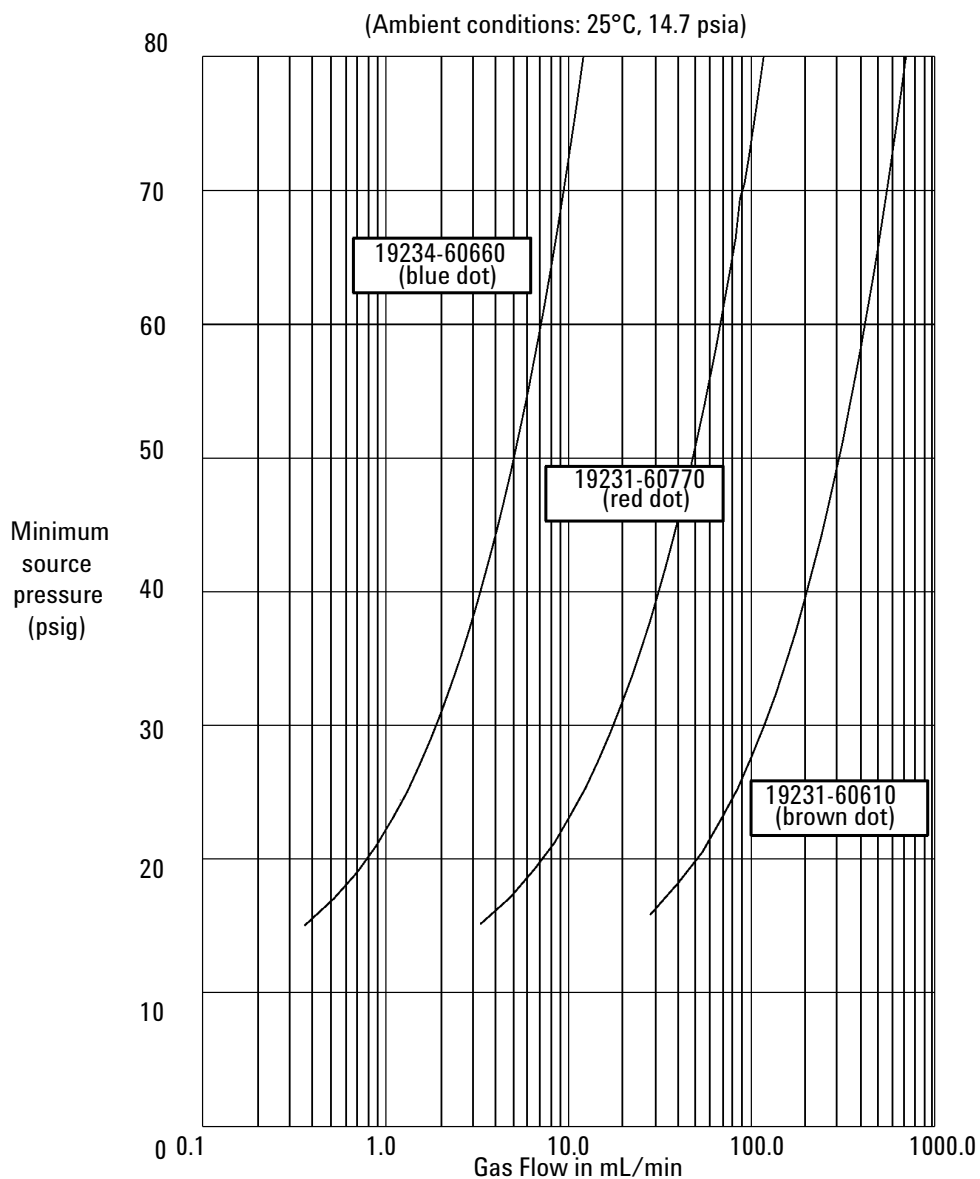


Figure 1. Pressure requirements for AUX EPC flow restrictors with air, nitrogen, or helium

Installing an Auxiliary Pressure Control Module
Install the new module

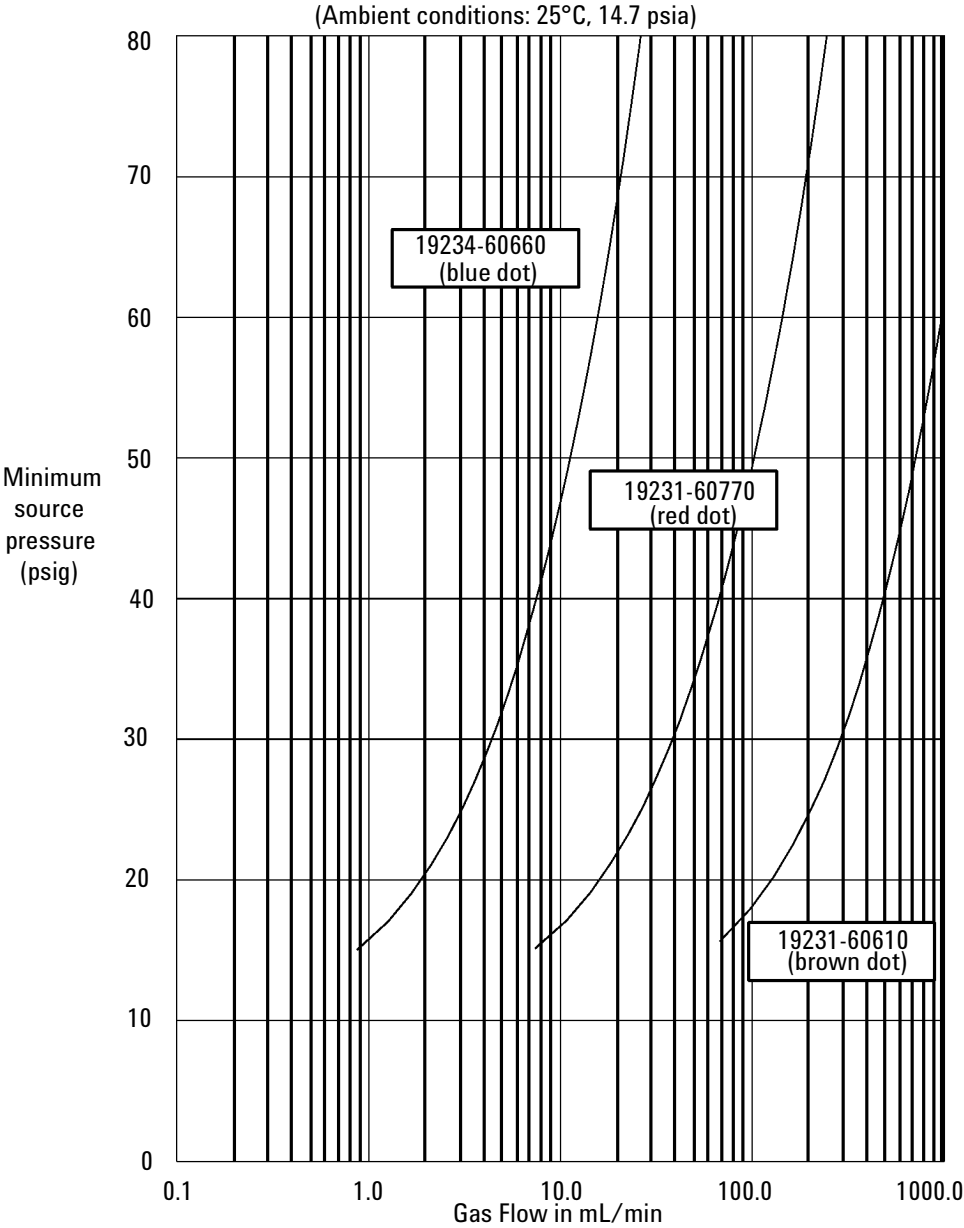
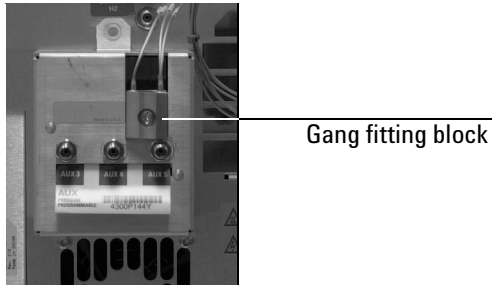


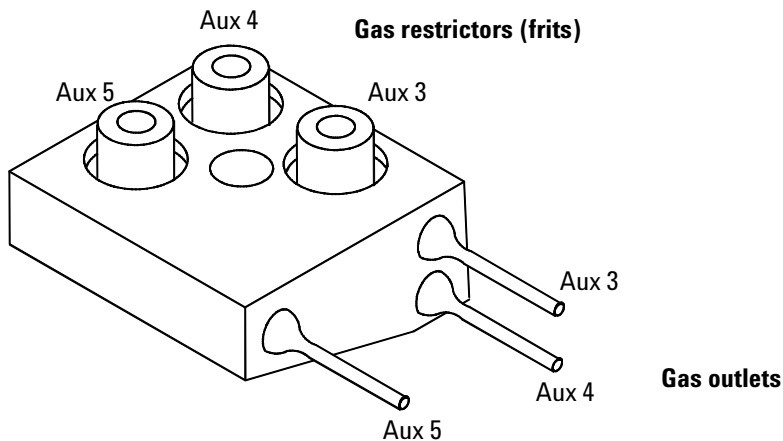
Figure 2. Pressure requirements for AUX EPC flow restrictors with hydrogen

Changing an auxiliary channel frit

1. Locate the gang fitting block. It connects the three gas outlet tubes for the auxiliary channels to the pneumatics module.



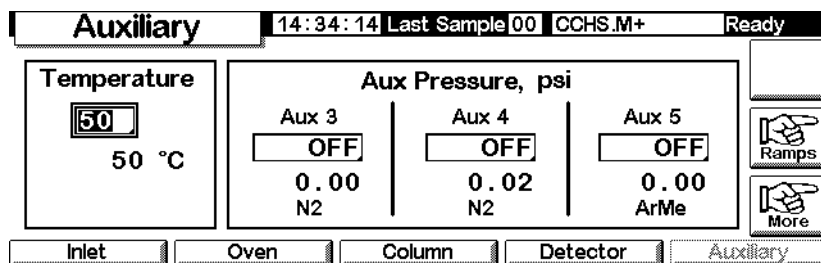
2. Remove the screw that holds the block to the pneumatics module. Pull the block free of the module and rotate it so that the frits are on top.



3. Pull the frit to be changed out of the block. Also remove the O-ring that seals it.
4. Place an O-ring on the new frit. Place the O-ring/frit combination in the block.
5. Reconnect the block to the pneumatics module. Tighten the screw firmly.

Restore the GC to operating condition

1. Reconnect all cables.
2. Restore power.
3. Display the Status / Settings / Auxiliary screen on the control module. If the module is installed correctly, this screen appears.



An actual flow value appears when gases are off or not connected. This is not an error. After the gases are connected and the detector is operational, the actual flow values are equal to the setpoint values.

If the screen reads Aux not installed, recheck your cable connections.

4. Enter 0 in each Aux field (3, 4, and 5) to zero the pressure sensors.
5. Connect the source gas lines to the module.
6. Restore carrier and other gases to the module and instrument.
7. Leak check all fittings.

Installing an Auxiliary Pressure Control Module
Restore the GC to operating condition



Agilent Technologies

© Agilent Technologies, Inc. 2004

All Rights Reserved. Reproduction, adaptation, or translation without permission is prohibited, except as allowed under the copyright laws.

Part number G3349-90007
First Edition, May 2004
Printed in USA

Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610

Acknowledgements

Torx is a U.S. registered trademark of Textron, Inc.

Safety Symbols

Warnings in the manual or on the instrument must be observed during all phases of operation, service, and repair of this instrument. Failure to comply with these precautions violates safety standards of design and the intended use of the instrument. Agilent Technologies assumes no liability for the customer's failure to comply with these requirements.

In the manual

A warning calls attention to a condition or possible situation that could cause injury to the user.

A caution calls attention to a condition or possible situation that could damage or destroy the product or the user's work.

On the instrument



See accompanying instructions for more information.



Indicates a hot surface.



Indicates hazardous voltages.



Indicates earth (ground) terminal.



Indicates explosion hazard.



Indicates radioactivity hazard.



Indicates electrostatic discharge hazard.



Indicates pinch hazard.



G3349-90007