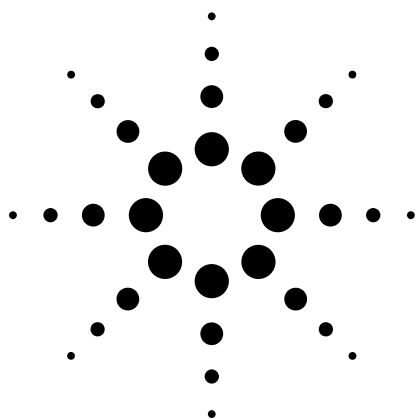




Installing the GPIB/RS-232 MIO Card into the 35900E Dual Channel Interface

The G1847A, GPIB/RS-232 MIO card allows the host computer to communicate with the 35900E through either an RS-232 or a GPIB connection.

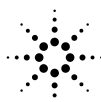
WARNING

To avoid personal injury, turn off the power and remove the power supply cord from the unit.

Caution

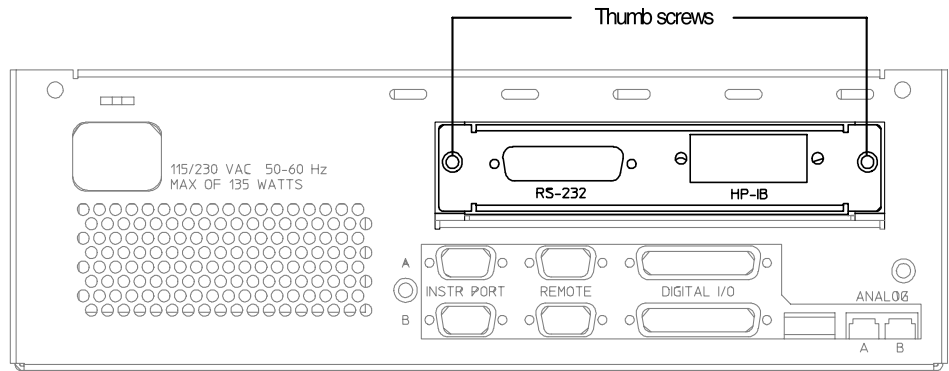
To avoid damaging this card during installation, use the following precautions:

- Use antistatic equipment during installation, such as a grounding mat and wrist strap.
- Hold the board by its edges.
- Do not touch the card's electrical components or edge connectors.



Agilent Technologies

Innovating the HP Way

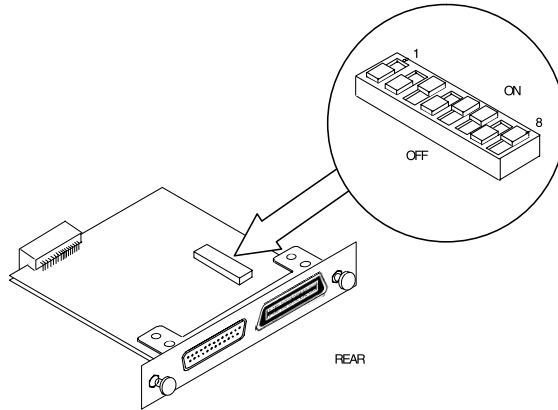


Steps

1. Turn off the power, and remove the power supply cord from the unit. Follow the appropriate configuration instructions that follow these steps.
2. Loosen the thumb screws and remove the installed card or cover on the rear of the 35900E.
3. Slide the new configured card into the unit. Be sure the edges of the card are in the guide rails.
4. Secure the card with the two thumb screws.

Configuring an GPIB/RS-232 MIO Card

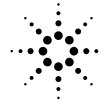
To configure the card for your system, you may need to change the settings of the switches located on the top of the card.



GPIB Configuration

When the position 1 switch is set to OFF, the MIO card uses the GPIB communication mode. In this mode, the ON switch position for switches 2 through 8 are defined in the following table:

Switch	ON Action
2	N/A
3	SRQ Enable
4	Add 16 to GPIB address
5	Add 8 to GPIB address
6	Add 4 to GPIB address
7	Add 2 to GPIB address
8	Add 1 to GPIB address



RS-232 Configuration

When the position 1 switch is set to ON, the MIO card uses the RS-232 communication mode. In this mode, the ON switch positions for switches 2 through 8 are defined in the following tables:

Switch	ON Action
2	Handshake Enabled
3	Hardware Handshake
4	Parity Enable
5	Even parity
6	Baud rate
7	Baud rate
8	Baud rate

Baud Rate	6	7	8
300	OFF	OFF	OFF
600	OFF	OFF	ON
1200	OFF	ON	OFF
2400	OFF	ON	ON
4800	ON	OFF	OFF
9600	ON	OFF	ON
19200	ON	ON	OFF
38400	ON	ON	ON

