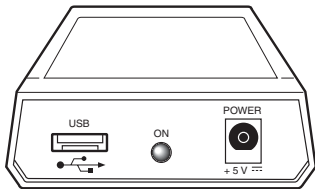


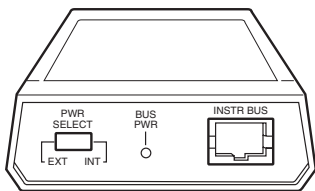
# BioLogic DuoFlow System USB Bitbus Communicator Catalog Number 760-2030

Dear Customer,

The BioLogic DuoFlow System USB Bitbus Communicator is the communication link between the DuoFlow Controller (computer), the DuoFlow Workstation and associated peripherals such as the DuoFlow Maximizer, Signal Import Module (SIM), Econo Gradient Pump (EGP), and fraction collectors.



FRONT VIEW



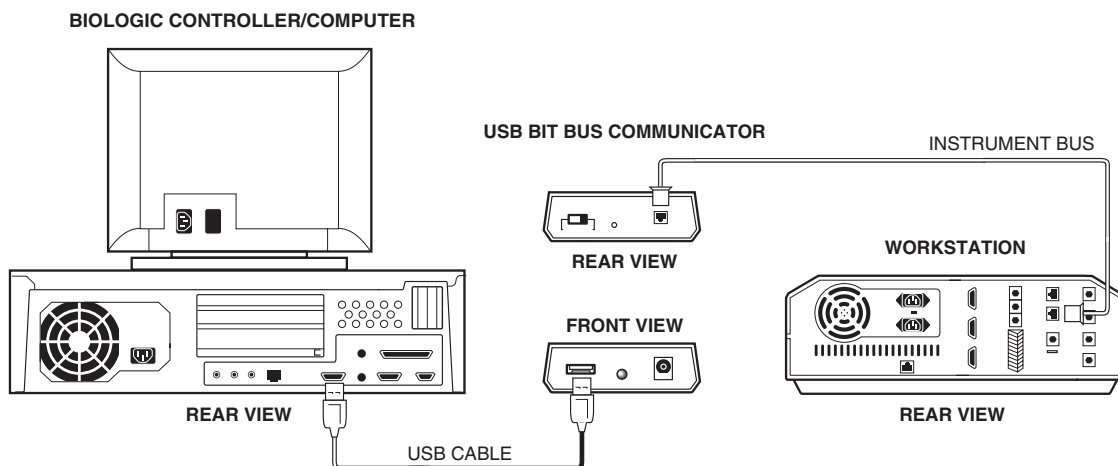
REAR VIEW

The USB Bitbus Communicator is powered directly from the Controller/ computer via the USB cable. In this case, the PWR Select switch should be set to INT.

If one or two SIM devices are used, the USB Bitbus Communicator must be powered by connecting a universal AC/DC in-line adaptor (Catalog number 760-2034). The PWR SELECT switch should then be set to EXT.

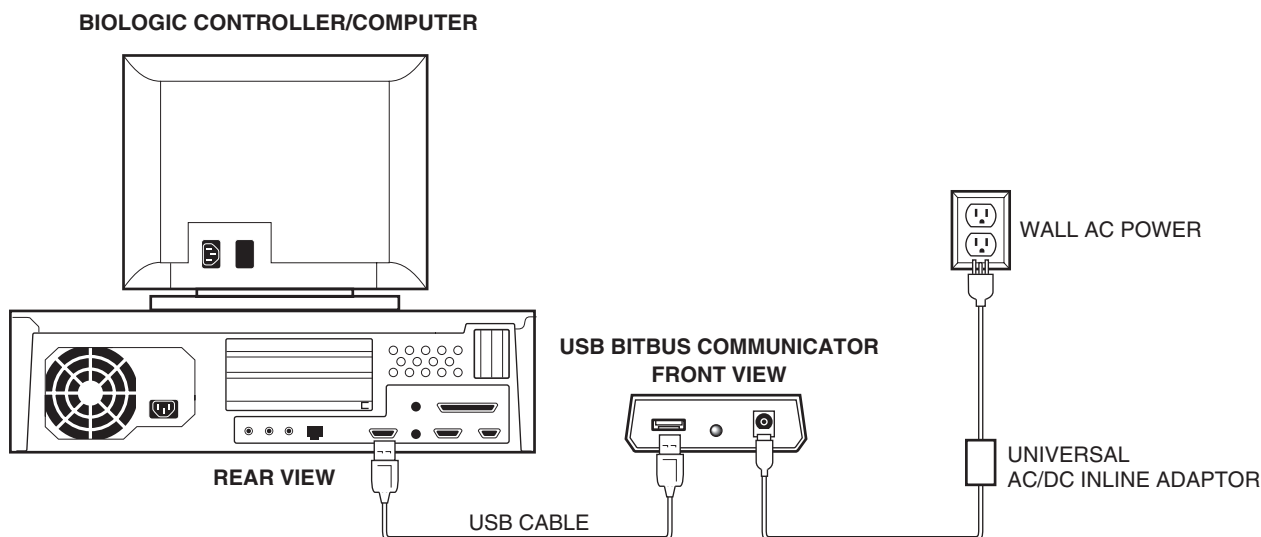
## Installation Instructions

1. With the power switched off, connect the USB cable between an available USB port on the Controller/ computer (front or rear) and the USB port on the USB Bitbus Communicator. (front view)
2. With the power switched off, connect the instrument bus cable between the Workstation and the INSTR BUS port on the USB Bitbus Communicator. (rear view)
3. The PWR Select switch should be set to INT.
4. Connect peripheral devices such as fraction collectors and auxiliary pumps to the Workstation via instrument bus cables or the Aux connector.

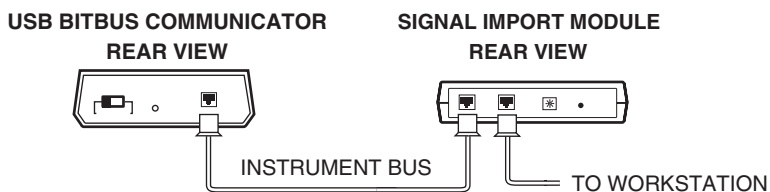


## Installation with SIM Devices

1. With the power switched off, connect the USB cable between an available USB port on the Controller/computer and the USB port on the USB Bitbus Communicator. (front view)
2. The PWR Select switch should be set to EXT.
3. Connect the universal AC/DC in-line adaptor (Catalog number 760-2034) to the USB Bitbus Communicator. When the AC/DC in-line adaptor is plugged into a wall socket, power is supplied to the USB Bitbus Communicator.
4. Switch on the Controller/computer and allow the Windows operating system to launch.
5. Verify that the BUS PWR light on the USB Bitbus Communicator is illuminated.
6. Connect one or two SIM devices to the USB Bitbus using the INSTR BUS port and cable. If two SIM devices are connected, ensure that one is set to device #1 and the other to device #0 using the small selector switch.
7. With the power switched off, connect the Workstation instrument bus port to a SIM device.
8. When all devices are connected, turn on power to the Workstation.



### 1 CONNECTING A SINGLE SIM



### 2 CONNECTING TWO SIMs

