



Connecting a *dataTaker* DT80 Range Data Logger to a Local Area Network (LAN)

1 Prerequisites

- Nil

2 Required Equipment

- *dataTaker* DT80 range data logger
- Ethernet cable
- Free network port on an active wall-plate, switch, router, etc.
- PC with free USB port and DeTransfer software (static IP address only)

3 Process

3.1 Physical Connection

1. Apply power to the *dataTaker*
2. Plug one end of the Ethernet cable into the dataTaker and the other end into a free, active Ethernet port (wall plate, router, switch or hub) so that the dataTaker is connected to the network. Ensure that both plugs 'click' into their sockets.



Figure 1 - Ethernet connection to the dataTaker

3. Now that the dataTaker is plugged into the network, it needs a network (IP) address.
 - a. If the network supports dynamic IP addressing it will automatically assign an IP address to the logger and you may proceed to the "Dynamic IP Addresses" section.
 - b. If the network does not support dynamic IP addressing then you must configure the IP settings of the dataTaker manually, proceed to the "Static IP Addresses" section.



3.2 Obtaining an IP Address

3.2.1 Dynamic IP Addresses

If the network has assigned the logger an IP address then it will be displayed on the “Eth IP” screen (on the dataTaker) as shown in Figure 2 (use the UP/DOWN arrows on the front of the logger to navigate to this screen).

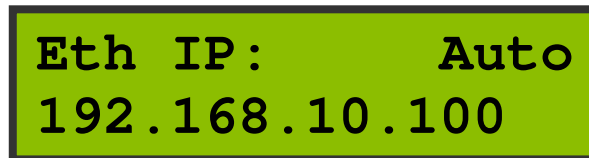


Figure 2 –‘Eth IP’ sample screen (Auto)

NOTES:

- If the IP address is displayed as 0.0.0.0, then the dataTaker is still looking for an IP address.
- If the IP address begins with ‘169.254’ then the logger has assigned itself an IP, not the server. In this case a static IP address must be used.

Once the logger has successfully been assigned an IP address, you will now be able to browse to the *dataTaker* web interface using any web-browser.

3.2.2 Static IP Addresses

To assign a static IP address you will need to follow these steps:

1. Obtain TCP/IP settings from the network administrator; you will need an IP address, subnet mask, gateway and DNS server addresses.
2. Plug the USB cable into the USB slave socket on the side of the dataTaker and the other end into a PC

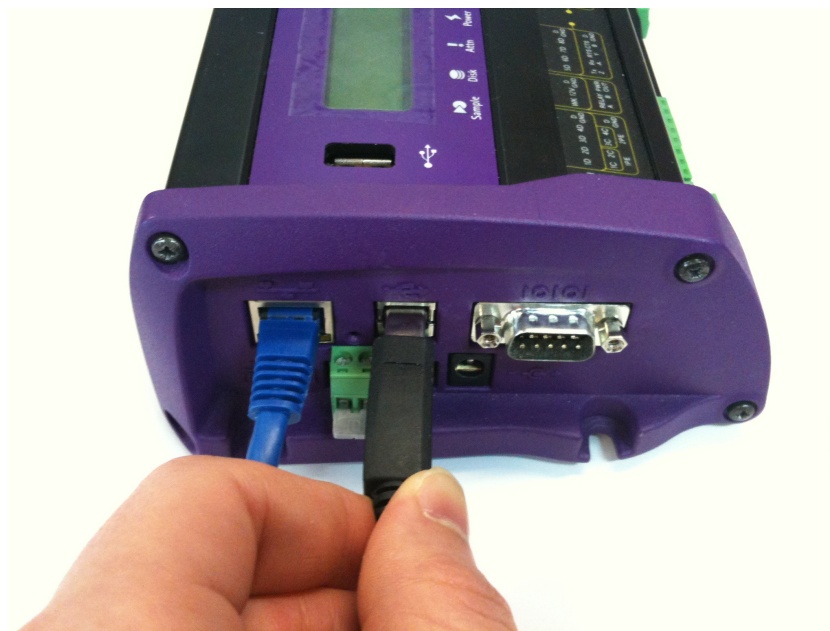


Figure 3 - USB slave connection to the dataTaker

3. Using DeTransfer, establish a connection to your dataTaker
4. Enter the following commands into the command window:



```
PROFILE"ETHERNET" "IP_ADDRESS"="XXX.XXX.XXX.XXX"  
PROFILE"ETHERNET" "SUBNET_MASK"="XXX.XXX.XXX.XXX"  
PROFILE"ETHERNET" "GATEWAY"="XXX.XXX.XXX.XXX"  
PROFILE"NETWORK" "DNS_SERVER_1"="XXX.XXX.XXX.XXX"  
PROFILE"NETWORK" "DNS_SERVER_2"="XXX.XXX.XXX.XXX"
```

(Where XXX.XXX.XXX.XXX are provided by the network administrator)

5. You will be able to check the IP address any time by looking at the “Eth IP” screen (on the dataTaker) as shown in Figure 2Figure 4 (use the UP/DOWN arrows to navigate to this screen).

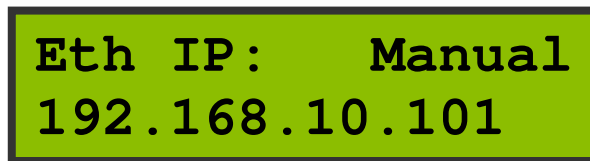


Figure 4 – ‘Eth IP’ sample screen (Manual)

Once you have successfully entered the TCP/IP settings, you will be able to browse to the dataTaker web interface by typing the address into a web-browser.