TN-0042

Securing the dataTaker DT80 Series Web Services

1 Settings changes (dEX)

NOTE: The changes specified in this document require your logger to be running firmware version 9.08 or above.

If your logger was configured using dEX, then the security settings are all located on security tab (click on the model number at the top of the tree, then Security tab).

Site information	Security
Regional Statistics CEM	To protect the logger from unauthorised access enter the credentials in the fields below. Additionally, any server that is not required can be disabled.
Power	Servers
Security	Command server 🗹 Port 7700
USB port	FTP server 🗹 Port 21
Sensor port	HTTP server ✓ Port 80
Ethernet	Modbus server ✓ Port 502
Email	Credentials for PPP and FTP
Modem	Username DATATAKER
Modem DDNS	Password ******** [] (mask)
Start up Keypad functions LCD	FTP Allow anonymous FTP ☑ Note: anonymous FTP is read only
Calibration	Command port Password (command password) ✓ (mask) Delay before locking 600 → seconds Note: when a password is defined, the logger automatically locks the command interface after this period of inactivity

1.1 Password protect the inbuilt FTP server

The FTP server is one of the two key methods for changing settings on your logger. The most important change when locking down the dataTaker web services is to prevent unauthorised access to this server. The recommended action here is to change the default FTP username and password and disable anonymous access.

1.2 Password protect the command port

Adding a password prevents your logger from accepting unauthorised commands. Once a password is set and the user is signed off the logger will not respond to commands. To access the command interface on a password protected logger, you must type the password then press enter. The command window is available within in the 'monitor the logger' interface.

The delay before locking period determines the timeout on the command interface, when this time expires the user will need to re-enter the password.

Page 1 of 3 TN-0042-A2



Technical Note

TN-0042

2 Settings changes (Command Language)

The below commands are equivalent to those mentioned in the previous section and should be put within a jobs code so that they are re-instated if the power to the logger is cycled

PROFILE FTP_SERVER ALLOW_ANONYMOUS=NO
PROFILE FTP_SERVER USER=DATATAKER
PROFILE FTP_SERVER PASSWORD=NEWPASSWORD
P14=120 'delay before locking
PASSWORD="MYPASSWORD"

NOTE: The above commands.

3 Limit access to the 'Monitor the Logger' interface

The *monitor the logger* interface is configured from the *Customise dEX* menu on logger home page.

3.1 Hide certain diagnostic features to keep the FTP password safe

Enter the settings as shown on the right for each of the menus. These will remove some parts of the web interface where a user might be able to view profile settings and see the FTP password.

The local documents link will not work because we have removed the FTP anonymous user so it should also be removed.

Save the changes by clicking Save Settings

3.2 Prevent editing of mimics

After you have added all of the mimics you wish to use in the web interface, you will want to prevent users from changing them. To do this, you will need to re-enter the *Customise dEX* menu and untick the box next to **Allow changes to mimics** in the *Security* page.

Save the changes by clicking Save Settings

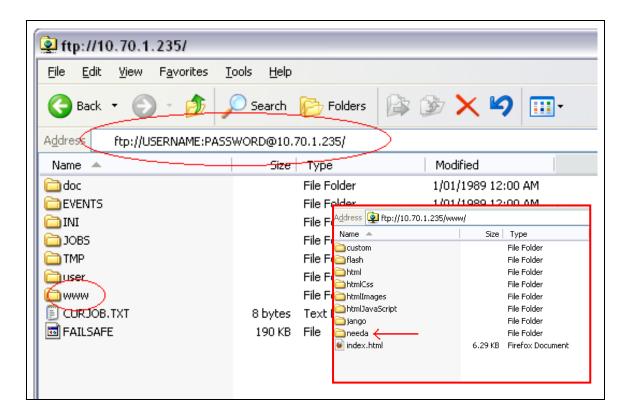
Status menu items
Include 'Welcome' menu 🗹
Include 'Schedule' menu ✓
Include 'Storage' menu 🗹
Include 'Diagnostics' menu
Include 'Program' menu
Include 'Parameters' menu
Include 'Profile' menu
Include 'Switches' menu
Retrieve menu items
Include 'Basic retrieve' menu 🗹
Include 'Advanced retrieve' menu ✓
Command menu item
Include 'Command' menu
Help menu items
Help menu items
Help menu items Include 'This interface' menu ✓
Help menu items Include 'This interface' menu ✓ Include 'Video tutorials' menu ✓
Help menu items Include 'This interface' menu Include 'Video tutorials' menu Include 'Local documents' menu □
Help menu items Include 'This interface' menu ✓ Include 'Video tutorials' menu ✓ Include 'Local documents' menu Include 'Datataker web-site' menu ✓
Help menu items Include 'This interface' menu ✓ Include 'Video tutorials' menu ✓ Include 'Local documents' menu ☐ Include 'Datataker web-site' menu ✓ Include 'Technical support' menu ✓
Help menu items Include 'This interface' menu ✓ Include 'Video tutorials' menu ✓ Include 'Local documents' menu ✓ Include 'Datataker web-site' menu ✓ Include 'Technical support' menu ✓ Lock mimics Allow changes to mimics □
Help menu items Include 'This interface' menu ✓ Include 'Video tutorials' menu ✓ Include 'Local documents' menu ✓ Include 'Datataker web-site' menu ✓ Include 'Technical support' menu ✓ Lock mimics

Page 2 of 3 TN-0042-A2



4 Hiding the 'Logger Home' menu items

Now that all of the settings have been saved, you will want to hide the *Customise dEX* menu. To do this, open Windows Explorer and log into your dataTakers FTP server. Use the IP address of your logger, user name and password you created in the profile earlier.



Once logged in, open the **WWW** folder and rename the folder 'needa' to anything else (eg. needa1234). Next time the logger home screen is shown the **Customize dEX** menu will not be shown and will not be accessible (it may be necessary to clear the browser cache for this page to reload correctly). It is also possible to hide the **Configure the logger** menu item using the same process but renaming the folder "jango" instead of "needa".