

DELOAD v2.37 RELEASE NOTES

Upgrade procedure

To reinstall DeLoad, first make sure it isn't running then uninstall using the windows Add/Remove Programs control panel. Previous settings for dataTakers and options are saved across to the new installation.

Installation procedure

You must be logged in as a user with administrative rights for nt, win2000 and win xp to install DeLoad
Shut down all programs including anti-virus software before installing.

1. Unzip the file DeLoad-2v37.zip to a temporary directory.
2. Run setup.exe

During installation, if you are prompted to keep existing files that are newer than those being installed by DeLoad, click yes to keep the existing newer files. If you keep getting a message saying that windows has to update files and you need to restart windows, start windows in safe mode and then install DeLoad.

You can have multiple installations of DeLoad. During the installation, just choose different installation directories for each separate installation. Each installation will track it's own set of dataTakers and settings.

The installation directory path is written to the windows registry whenever deloadop.exe is run and if there is no install.ini file present in the directory in which deloadop.exe was run from. DeLoad reads this path every time it starts so that running DeLoad from a batch file still works. If this path becomes incorrect because you ran deloadop.exe inadvertently from a location other than the installation directory, (ie: you copied deloadop.exe somewhere else), (delete the newly created install.ini from the wrong directory and re-run deloadop.exe from the correct installation directory), twice, to re-create the correct registry entry.

Known issues

ISSUE: If a modem connection is lost during an unload a NO CARRIER message will appear at the end of the data.

WORK-AROUND: For some modems you can put an ATQ1 in the modem init string. Beware that we have found that for some PSTN modems this has caused problems such as not being able to dial more than once, so you must try this with the particular modems first.

Software Change History

Version 2.37, 05 May 2008 (this release)

=====

- FIX: Fixed a bug where the decimal point separator for the ulast timestamp saved to disk reflected that set in the regional settings for windows. This caused all data to be unloaded when ULAST was selected for unloading. dataTakers can only deal with the decimal point separator, and so DeLoad had been changed so this is always the case despite the regional settings.

- FIX: This version fixes having to install DeLoad in safe mode when windows keeps asking to restart the systems so that files can be updated. The following files have been removed from the installation.

olepro32.dll
stdole32.tlb
oleaut32.dll

These files are part of windows default installations, however if this version won't install use version 2.36. DeLoad versions 2.36 and 2.37 beta are identical in all other respects.

Version 2.36, 25 April 2007

=====

- ADD: Added support for DT85 dataTakers.

Version 2.35, 18 April 2007

=====

- ADD: DeLoad now creates a file, one for each dataTaker schedule, named currentfile.txt, where n is the schedule letter, in the directory where dataTaker data is being unloaded to. These files contain the file name of the file that data is currently being unloaded to. The files are updated each time data is unloaded from a dataTaker using DeLoad.

Version 2.34, 20 September 2006

=====

- FIX: Data files are now opened or created after the first crlf is received from the logger rather than just before the unload command is sent. Before, empty files could be created if the connection to the logger dropped out between opening a file and data being returned.

- FIX: Fixed a bug where unload from all dataTakers returned the connection name from found.ini rather than from the dataTakers found by autobaud.

- FIX: Data file time stamps are now in 24 hour format. Before they were in 12 hour format meaning that am files could get overwritten if the pm time was the same.

- FIX: Append to file now works in all situations, including saving data to mdbs and for dlr files. When saving to mdbs, data is stored to temp files in unloadroot as well as to the schedn files. Then when data is added to an mdb, the temp files are used which contain just the data for the current unload.

- FIX: Checked multiple installations - this actually all works but the install.ini file in the shipped distributions had text in it. DeLoad relies on this file being empty the first time it's run so that the appropriate installation root registry setting is written. Then after that each installation of DeLoad looks for a registry setting named by the current directory it is run from, and makes that its working directory.

- CHANGE: For modem connections, DCD is now checked to make sure it's inactive before an attempt is made to connect to the local modem. If not, an error occurs.

- CHANGE: An error message is returned if a modem connection drops out during an unload, and unloading is canceled. Before, you had to wait the connection timeout.

- CHANGE: Modem connections now use DCD to determine if a remote modem is connected to.

- CHANGE: DeLoad now has a log file. You can turn it on and off and set the max size in Options. When the max size is reached, the log file is backed up once. The next time it's overwritten. A button has been added to options to view the log file.

Version 2.33.1, 18 August 2006

=====

- FIX: Fixed a bug where for modem connections the ATH hang up command was not being sent. This hasn't seemed to be an issue in most cases because when the comport is closed most modems automatically drop the connection. Problem known to affect 905U-D radio modems.

- FIX: Fixed a problem whereby DeLoad would appear to hang when trying to download data with missing characters (eg: over a bad modem connection). Caused by a 4 second timeout for each missing character when converting a DT80/800 series file to csv. Changed the way the the csv conversion process works so as to remove the timeout.

- FIX: Added a check on the date/time stamp for records being converted to csv format for DT80/800 series. Records with a bad date/time are not copied to the csv file.

-FIX: Added a check on the record length for records being copied to csv format for DT80/800 dataTakers. If the actual length doesn't agree with the length at the end of the record, the record isn't copied to the csv file.

- FIX: Added a check on the data/time stamp written to file for ulast unloads for DT80/800 series dataTakers. If the date/time stamp is of the incorrect format it isn't written to file. When read from file, if it isn't the correct format the next unload will unload all data with no from data/time stamp.

- FIX: Fixed the last missing character problem for the data time stamp saved to file – now the time will always have 8 characters for the seconds.
- FIX: You can now have commas in the phone number and initialisation fields for modem connections. Commas are saved to file as | s and converted back to commas when the file is read.
- FIX: Fixed a bug where the connection timeout was nothing after you did a find dataTakers. Now if the connection timeout field is ever blank it is automatically set to 10.
- CHANGE: Changed the progress bar so that it only updates for good records that are converted to csv format. This gives a visual indication of how many bad records have been removed.
- CHANGE: Changed the delay between sending a +++ and commands after that for modems from 100 mSecs to 1.5 seconds.

Version 2.30, 10 July 2006

- FIX: Fixed installer bug introduced in version 2.29 which resulted in DeLoad not being able to be installed. Caused by missing files in the installation.
- FIX: Fixed a bug for DT500 series dataTaker CSV unloads where data records were being doubled up.
- CHANGE: Moved the ConnectionTimeout parameter from the options.ini file to the found.ini file and added a text box in the SetCom form allowing this parameter to be individually set for each dataTaker from DeLoad Options.
- CHANGE: Removed check for BEGIN"JOBNAME" at start of programs being sent to DT80/800 series dataTakers. Now single line programs with out BEGIN and END statements can be sent.
- CHANGE: If the 'Show Unload Form' check box in DeLoad Options is unchecked, forced 'Unload data when DeLoad Starts' Check box to be checked. Otherwise DeLoad starts and sits in the background doing nothing when both boxes were unchecked.
- CHANGE: Removed the 'Hide DeLoad after unload' check box in Unload Options because it actually didn't do anything.

Version 2.29

- FIX: Removed the time stamp at the start of dlr lines of data for DT500 Series dataTakers to match the change in version 2 of Deview.

- ADD: Added support for DT81s.

Version 2.27

- FIX: Fixed up some problems with sending emails where the wrong data files were being sent.
- FIX: Made save to one file work properly with files sent via FTP.

Version 2.26

- FIX: deload 2.25 did not work properly for dt80/800s
- FIX: fix stretched panel on datataker options form
- FIX: fix bug in del job function for dt80/800s
- CHANGE: change the way job lists for dt80/800s were gotten to cater for differences in version 4 and 5 of dt80/800 firmware - now uses dirjob* to get job details list
- CHANGE: changed the file extension for unloaded non csv data files where append to last saved file is ticked - now it's .txt and the dlr header isn't included - will fix this later - means that these files are just text data in format mode and won't work with deview or delogger

Version 2.25

- FIX: Com ports over 8 didn't seem to work properly in some circumstances so this was revised and fixed.
- CHANGE: Changed the progress bar for DT800s to count records rather than characters using dirjob* rather than dirtree.
- CHANGE: Removed the 'check that Q was received' check before unloading data and put 5 blind Q sends without a check.
- ADD: Added DT80 support.

Version 2.24

- CHANGE: Added a check when saving the ULAST date/time stamp to file that only writes the file if the length of the data/time stamp is correct. This is because if data returned by a logger is corrupted then if the data/time stamp is bad the dataTaker will

never unload using ULAST again as the U command fails and there is never any new data to get the next date/time stamp from. The actual contents of the date/time stamp is not checked as this means knowing the date and time separator characters meaning the P values have to be gotten etc, - too much mucking around. An incorrect length will catch corruption in 99.9 % of cases probably.

Version 2.23

=====

- FIX: Fixed save to dlr format by including [DT500 timestamp + difference bewtween dt basedate and ms basedate] for DT500 dataTakers
- FIX: Fixed save to dlr format by modifying records with 012345 and working out the new checksum for DT800 dataTakers.
- FIX: Fixed missing header information by adding S,012345 to the start of the header line. Also, changed the header from STATUS10 to STATUS14 information, and included the new checksum.
- FIX: Ulast for non csv data or dlr format data can now be performed. This inherently fixed a slight bug where for DT500 datatakers, if there was only one record after the ulast timedate stamp, that record wouldn't unload until the next unload.

Version 2.22

=====

- FIX: datatakers with CEMs never get recognised becuase of a slight bug when deciphering the results of TEST0
- CHANGE: made the connection timeout applicable to all parts of modem connections - before, sending TEST0 to check for a logger used delay modem connections when using the terminal in deload options were always set to 57600 instead of what the user chooses in the setcom form - this was stupid and is now fixed.

Version 2.21

=====

- FIX: fixed a bug where finding a memory card using a DT500 didn't work when erasing the card
- FIX: added delogger dlr headers to non-csv data files
- CHANGE: removed dtx terminal from the installation. datataker.ocx terminal will be added later

Version 2.20

=====
- FIX: fixed a bug where if autobaud worked but the datataker id couldn't be determined (if a password was given to datataker) then if the enter password box was cancelled the progress bar form didn't come back and a non-existent logger was found

- CHANGE: also changed the delay when waiting for the datataker id in rs23autobaud to wait 1 sec longer

Version 2.19

=====

- FIX: fixed a bug where if com 1 wasn't the first port then finding all dataTakers didn't work

Version 2.18

=====

- CHANGE: changed deload.exe so that it gets the current switch settings via STATUS9 and P22 and sets these back to how they were before unloading data using deload

- CHANGE: changed the email subject so that a list of schedule descriptions is appended after the site id description when all schedules isn't used - for all scheules just the site is put in the subject

Version 2.17

=====

- ADD: added a simple terminal program as a separate exe with a start menu icon

- ADD: added a dialog box for creating an icon which allows the icon name and icon function to be set

Version 2.15

=====

- FIX: fixed a bug where dragging a file onto an icon made getfilesize not work coz of the quotes in commandlineargument causing the progress bar to not work properly.

- FIX: fixed a bug where ftp was broken when the name of the ftp directory in which html files are kept was changed to description

- CHANGE: changed sending a program to ram and flash so that the file open dialog box opens before connecting
add the send command to datataker #commmand argument for delpoad.exe including the option to run in terminal windows #command#

- CHANGE: improved dt500 unloading by eliminating the doubled up record at the start of unloaded data caused by the ulast date/time stamp not being correctly increased by 1 second

Version 2.14

=====

- FIX: fixed a bug where if connecting to a modem timed out when sending a program to a logger, options form was still running invisible

- FIX: fixed a bug where deload.exe would probably only work with datatakers on comports under 9

- FIX: fixed a bug where when downloading from a datataker over a slow modem link the results of dirtree to get schedule data sizes didn't work properly - caused by a vb timing issue - fixed by clearing the character window after inbetween doing a dirtree a: and dirtree b:

- FIX: ulast unloading of data was doubling up the first record between unloads for dt800s - fixed this by checking the time and date for every record (converted to csv for dt800s only) against the lastunloadeddetails time and date and dropping the record if it already has been downloaded

- FIX: the first unload "with save data to one" file ticked only downloaded the header the first time - this was fixed - caused by an err not being cleared

- FIX: fixed a bug where opening the setcom form for a dataTaker made an existing desktop icon not work any more

- CHANGE: deload options now finds com ports automatically when started and when the find button is clicked

- CHANGE: increased the number of rs232 com ports from 8 to 20 (this should have been easy but it was really hard because found loggers were saved as eg: 13458 :())

- CHANGE: added a windows registry entry which contains the installation directory for deload - the current directory is now looked up from here rather than using CURDIR meaning that if deload is run from a batch file in another directory everything still works

- CHANGE: added a modemtimeout parameter in options.ini - this because mobile phone modems take longer to connect

- ADD: added dt53 to deload

Version 2.13

=====

- FIX: fixed a bug where if deload.exe was clicked (no arguments) and there were multiple dataTakers, the first dataTaker that wasn't found caused deload to exit thereby not processing the rest of the dataTakers - this was a stupid bug if comandlinearguement <> ""

then exit sub was actually if comandlinearguement = "" then exit sub

- CHANGE: change the data type for data fields in mdbs from string to double

Version 2.12

- FIX: the state of the mdb tick box was not being saved and loaded properly resulting in the user settings being changed to no tick - this also resulted in csv being unticked when mdb was ticked causing the copy data to database function to crash - mdb tick boxed state fixed and check put at start of copy to database procedure that exits if csv for that schedule is not true

- FIX: when a datataker was deleted the email list file wasn't because the datataker tag used was the number in the drop down list rather than the actual tag ie ::45

- FIX: added an extra check for no data for dt800s where [end of schedule, end of unload] and [H line, end of schedule, end of unload] are all catered for - before it was possible that if a halt hadn't been sent deload could think that there was data when there wasn't in particular when there was only one line of data

- FIX: removed shift A to send all text in the terminal transmit window - capital A couldn't be typed in :(

Version 2.11

- FIX: changed sending programs to flash so that lines starting with a '\', 'D='\ and 'T='\ are not sent and added the # character to allowed characters (it was left out meaning eg : 1#L wouldnot be sent), and so that a RESET can be sent if it isn't at the start of the line eg : K1="RESET"

Version 2.10

- FIX: Fixed a bug when sending a program to flash where lines with a RESET in them aren't sent and lines starting with a tab aren't sent.

- ADD: DeLoad has been changed so that now it checks DT800 records as they are converted to CSV format for correct record length. If the length is incorrect the record is dropped and won't appear in the CSV file. If any bad records are found a file is created in the unload directory called LostDT800CSV.txt, in which the time, date, DT800 description and total number of records lost for all schedules for the last unload for that dataTaker is written. This check is only done if CSV is ticked and is only for dt800s.

Version 2.09

- FIX: Fixed a bug where html data would be doubled up if the number of records unloaded since the last unload was less than the record window.
- FIX: Discovered that ftp doesn't work with the Godzilla ftp server. This issue wasn't resolved. Note that ftp has only been tested properly on a unix style ftp server such as the one that comes with freebsd unix.
- CHANGE: Modem connections have been improved by allowing the local modem baud rate to be chosen. This is for modem configurations where the baud rate of the local modem flows through to the remote dataTaker.

Version 2.08

- FIX: Fix a bug where email addresses and schedules didn't match properly.
- ADD: Added a "run datataker command after unload" function.
- ADD: The connection timeout can now be manually set.

Version 2.07

- FIX: Fixed a few bugs with the terminal windows where the transmit window didn't behaved as it should.
- ADD: Added the ability to specify individual email addresses for each datataker schedule. Also now you can add some text in the email body.
- ADD: Also added shift A to send all lines in the transmit window.

Version 2.06

- FIX: ulast didn't work properly if data was unloaded using deload whilst a datataker was logging, as the date,time stamp was written when the unload was started rather than taken from the last line of unloaded data. This has been fixed - now the ulast date,time is taken from the last line of data that was actually unloaded.
- FIX: End of unload detection has been improved such that no data is lost if connection to a dataTaker fails whilst unloading using Deload's ulast.
- FIX: Also fixed a bug in deloadop.exe where if you dragged a file onto an icon with an id for a datataker that doesn't exist, deload options would keep running with no form visible.
- ADD: Added a "run program after unload" function.
- ADD: A text based terminal window has been added allowing dataTaker text commands to be sent from Deload

Version 2.05

- ADD: Added the ability to type in a name for tables in the access database. Default is Schedulex.

Version 2.04

- FIX: connecting to a dataTaker from the options control panel when trying to send a file etc. did not recognise a dt51 with display (and probably a dt50 as well). This was a silly bug and was fixed.

Version 2.03

- FIX: got rid of the zero data record that gets appended to unloaded data when a H is sent

- FIX: for ulast, if the dt800 is logging data during the second that the time stamp is saved to file, then data from the point during the second that the timestamp was saved until the end of the second won't be unloaded next time - this is because ulast only works with increments of 1 second, not subdivisions of a second - the other way to do this is to unload all data for the second but this would result in doubled up data for the entire second

- ADD: added copy to mdb file tick box - only creates access 2000 database - access 95 and 97 to be added if requested by customers (one database per dataTaker, 1 table per schedule, 1 field per measurement type)

Version 2.02

- FIX: added a /m after u for dt500's because if logger was returning errors data unload details didn't work for each schedule and deload would not hang but stop when getting data unload details for sched a

- FIX: removed frmtaskbar.mscomm1.rthreshold = 0 just before status10 is sent to get headers - this was a stupid bug

- FIX: passwords didn't work at all - passwords were supposed to be saved in found.ini for each datataker but they ended up being saved in options.ini which is all wrong - changed things back to saving passwords in found.ini

- FIX: fixed up what happens when there is a password and when the one typed in is wrong ie : unknown datataker type etc.

- FIX: fixed a strange bug where when passwords are used a 2nd instance of deload.exe was started causing both instances to crash - didn't work

out what was going on but adding a msgbox just before the error occurred
fixed everything - hmm strange - this bug caused deload to crash
if there was a datataker password and a file was dragged onto a datataker icon

Version 2.01

- FIX: loading doesn't detect the absence of dataTakers
- FIX: auto click the apply button when unloading data from the datataker's unload options panel
- FIX: run time error on smtp test
- FIX: no help for save to disk text box
- FIX: help text for unload button for loggers is wrong
- FIX: help for ulast is wrong
- FIX: schedules frame help text is wrong
- FIX: edit path button help is wrong
- FIX: load buttons help text say a instead of the
- FIX: datataker list cbx help text is wrong
- FIX: testing dialling a datataker when no phone line is plugged in is all wrong
- CHANGE: change the name of smart.exe and options.exe to load.exe and loadoptions.exe
- CHANGE: memory card found message when loading a program - needs wait 1000 after it

Authorised distributor

In Australia:

For customer service, call 1300-735-292
To email an order, ordersau@thermofisher.com
To order online: thermofisher.com

In New Zealand:

For customer service, call 0800-933-966
To email an order, ordersnz@thermofisher.com
To order online: thermofisher.com

ThermoFisher
S C I E N T I F I C