

## DTUSB v1.02 RELEASE NOTES

---

---

Windows driver to provide TCP/IP access to DT80 Series logger over USB.

Requires Windows XP or later

### Installation

---

1. Uninstall any existing version of DtUsb, using the start menu link
2. Run DtUsbPackage32-102.exe (32 bit OS) or DtUsbPackage64-102.exe (64 bit OS)

### Contact Details

---

Thermo Fisher Scientific Australia Pty Ltd  
5 Caribbean Drive (PO Box 9092) Scoresby VIC 3179 Australia  
Phone: +613 9757 4477 Fax: +613 9753 3483  
Email: support@datataker.com.au

### Change History

---

---

DtUsb version 1.02 02-Jul-2010

---

DtUsb:

- no changes

UsbSerialiser:

- FIX: crash following serialisation on Windows 7 (#2333)

Installer:

- FIX: now includes multi-lingual version of DPINST driver installer, to allow installation on non-English Windows (#2337)

DtUsb version 1.00 16-Mar-2010

---

Initial release

## Authorised distributor

---

### **In Australia:**

For customer service, call 1300-735-292

To email an order, [ordersau@thermofisher.com](mailto:ordersau@thermofisher.com)

To order online: [thermofisher.com](http://thermofisher.com)

### **In New Zealand:**

For customer service, call 0800-933-966

To email an order, [ordersnz@thermofisher.com](mailto:ordersnz@thermofisher.com)

To order online: [thermofisher.com](http://thermofisher.com)

**ThermoFisher**  
S C I E N T I F I C