



# Plexor® Analysis Software, Forensic Release

---

Prior to analyzing data from the Plexor® HY System, the Plexor® Analysis Software (Forensic Release) must be downloaded and installed on a computer with Windows® 98 or newer operating system. This software is available free of charge and can be loaded on any compatible computer.

View the [system requirements](#) and a [version history](#) of the Plexor® Analysis Software

Upon completing the request for download, an email will be sent to you with a link to the downloadable files.

To download the Plexor® Analysis Software (Forensic Release), please complete the form below:

**Please provide your contact information:**

Underlined entries are required.

First Name

Institution

Street Address

PO Box

City

Country

Telephone

Email

Last Name

Department

Bldg/Room

Address is for:  Work  Home

State/Region

Postal Code

Fax

**If you recently moved, help us stay current - please tell us where you were before!**

Previous Institution

I give Promega or an authorized Promega distributor permission to contact me at the addresses that I provide.

Moved here

Select One

Submit