

## Licensing Information

This product includes copyrighted third-party software licensed under the terms of the GNU General Public License (Version GNU 2.0. The full terms of conditions you'll find at <http://www.gnu.org/licenses/old-licenses/gpl-2.0>).

Specifically, the following parts of this product are subject to the GNU GPL:

- The Linux operating system kernel
- The iptables packet filter and NAT software
- The busybox swiss army knife of embedded linux
- The zebra routing daemon implementation
- The udhcpd DHCP client/server implementation
- The pptp-linux PPTP client implementation
- The bridge-utils package
- The Ez-ipupdate dynamic DNS implementation
- this is just an incomplete example list

All listed software packages are copyright by their respective authors.

Please see the source code for detailed information.

## Availability of source code

Sartorius Lab Instruments GmbH & Co.KG. has created an Image file with the full source code of the GPL licensed software, including any scripts to control compilation and installation of the object code.

You can download that image from the sartorius web page free of charge.

Link to the download file: <https://www.sartorius.com/en/products/weighing/laboratory-balances/premium-level-balances/cubis/cubis-individual/q-app-center...../labpro-cubis-gpl.tar.gz>

All future firmware updates will also be accompanied with their respective source code.