

Magnis NGS Prep System Protocol Download Instructions

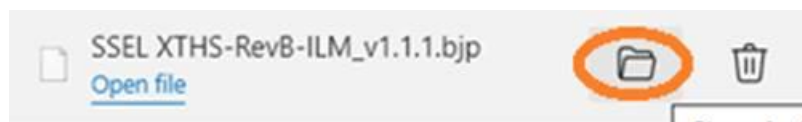
You can either download the protocol on USB and then upload to the instrument or download from Magnis touchscreen if connected to internet

Download the protocol on USB and then upload to the instrument

Tip! Use a USB stick that is unencrypted and FAT32 compatible. The USB stick should be blank, with no other saved files. Any files on the USB may be erased once inserted into the Magnis system.

Don't have the right USB stick?
Contact your local support team for assistance.

- Select the appropriate protocol from the Protocol Menu.
- Click **Download**, and once you see the protocol file, **Press** the “Folder” icon (see icon circled in orange below. Don't press “Open File”), save the file (*.bjp) to a USB stick



- Insert the USB stick with the saved file into an available port on the front of the Magnis
- From the home screen, press **Settings** > Protocol Update > Update Protocol for a protocol
- Select USB disk and select the file that you want to upload.
- Press **Import** for protocol
- Select **Update Protocol** from the **Protocol Update** screen.
- At the top of the **Protocol Update** screen, select **USB**.
- Press the USB Disk folder to expand it until you see the protocol file (*.bjp) downloaded from the webpage.
- Press the **protocol file (*.bjp)** to select it.
- Press **Import**. A message box opens notifying you that the protocol was imported.
- Press **OK** to close the message box. You are returned to the **Protocol Update** screen.
- Press **Close**. You are returned to the Protocols screen. The name and version number of the imported protocol is now listed on the screen.

Download protocols using the Magnis touchscreen

For Magnis systems with a network connection, new protocols can be downloaded using the unit's touchscreen interface.

- From the home screen, press Settings > Protocol Update > Update Protocol for a protocol
- Select Online Download and check the file that you want to download.
 - You may download more than one file at a time.
 - Download time may vary, depending on the connectivity speed.
- Press Download.
- For new protocols, you can return to the Protocols screen after the download completes to confirm that the new protocol is now listed.