

PLAT Recommended PC Setup

PC setup

CAUTION

Wrong settings

The recommended PC to use comes bundled with the Parallel Capillary Electrophoresis Instruments.

If the computer settings do not match that below communication issues to the instrument are possible, resulting in lost time and productivity.

- ✓ If a different PC is to be used or if any changes are made to the existing PC, check the following settings and enabled/disabled the PC settings to match the recommended setup.

Date and Time Settings

- 1 Go to **Settings> Time & Language> Date & Time** and set to Local Date & Time.
- 2 Go to **Settings> Time & Language> Date & Time** and turn off the option **Adjust for daylight saving time automatically**.
- 3 Win 10 - Go to **Settings> Time & Language> Region** and set the **Regional format** to **English (US)**.

Win 11 - Go to **Settings> Time & Language> Language & region** and set the **Regional format** to **English (US)**.

USB Power Management Settings

- 1 Go to **Device Manager > Universal Serial Bus Controllers > USB Root Hub (USB 3.0)**.
- 2 Right-click **Properties > Power Management** and deselect the **Allow the computer to turn off this device to save power** check box.

NOTE

To ensure the best operating conditions, it is recommended to close all other applications while the data acquisition software is running.

[System Overview](#)
PC setup

USB Power Management Settings

- 1 Go to **Device Manager > Universal Serial Bus Controllers > Intel® USB 3.1 eXtensible Host Controller (Win 10) or Intel® USB 3.2 eXtensible Host Controller (Win 11)**.
- 2 Right-click **Properties > Power Management** and deselect the **Allow the computer to turn off this device to save power** check box.
- 3 Win 10 - Go to **Settings > System > Power & Sleep** and select **Never**.
Win 11 - Go to **Settings > System > Power > Screen, sleep, & hibernate timeouts** and set all to **Never**.
- 4 Win 10 - Go to **Additional Power Setting** and select **High Performance**.
Win 11 - Go to **Power Mode** and select **Best Performance**.
- 5 Go to **Settings > System > Notification** and set all to **Off**.
- 6 Go to **Settings > Gaming** and turn off the Xbox Game Bar:
 - a Win 10 - Deselect **Open Xbox Game Bar**.
Win 11 – **Game Bar** set to **Off**.
 - b Set **Game Mode** to **Off**.
- 7 Win 10 - Go to **Settings > Privacy** and set all to **Off**.
Win 11 - Go to **Settings > Privacy & security > Recommendations & offers** and set all to **Off**.
- 8 Win 10 - Go to **Settings > Update & Security > Windows Update > Advanced options** and set all to **Off**.
Win 11 - Go to **Settings > Windows Update > Advanced options** and set all to **Off**.
- 9 Win 10 - Go to **Settings > Update & Security > Windows Security > Virus & Threat Protection > Virus & Threat Protection Settings** and set all to **Off**.
Win 11 - Go to **Settings > Privacy & security > Windows Security > Virus & Threat Protection > Virus & Threat Protection Settings** and set all to **Off**.
- 10 Win 10 - Go to **Settings > Update & Security > Windows Update** and select **Check Now/Refresh** to update to the latest available version.
Win 11 - Go to **Settings > Windows Update** and select **Check Now/Refresh** to update to the latest available version.
- 11 Win 10 - Go to **Settings > Update & Security > Windows Security > Core isolation > Core isolation details** and set **Memory integrity** to **Off**.
Win 11 - Go to **Settings > Privacy & security > Windows Security > Core isolation > Core isolation details** and set **Memory integrity** to **Off**.