



Configure Windows 10 Version 1703

- [MUST] 1. **System** (Microsoft Control Panel): Register Windows with Microsoft.
- [MUST] 2. **File Explorer Options** (Microsoft Control Panel): In the **View** tab:
- Select **Always show menus**.
 - Select **Display the full path in the title bar**.
 - Clear **Hide extensions for known file types**.
 - Clear **Use Sharing Wizard**.
- [MUST] 3. **Start > Settings > Update & security**:
Click **Check for updates** to check for updates and apply all critical security patches. Before proceeding, ensure that all updates are downloaded and installed. Ensure that there is no reboot pending.
- [MUST] 4. **Indexing Options** (Start > Windows System > Control Panel): Disable indexing.
Click the **Modify** button. Clear all drives and locations.
- [MUST] 5. **Start > Settings > System > Tablet Mode**: For **When I sign in**, select **Use desktop mode**.
- [MUST] 6. **Power Options** (Start > Windows System > Control Panel):
- a. As preferred plan select **High performance**
 - b. Click **Change Plan settings**
 - c. Set the option **Put the computer to sleep** to **Never**
 - d. Click **Change advanced power settings**.
 - e. Open the nodes for **Hard disk > Turn off hard disk after**.
 - f. Set the Minutes to 0 (=Never).
- [MUST] 7. **Start > Settings > Apps > Offline Maps**: Turn **Metered connections** and **Map updates** off.
- [MUST] 8. **Administrative Tools** (Start > Windows System > Control Panel): Configure security options.
- a. Double-click **Local Security Policy**.
 - b. Navigate to **Security Settings > Local Policies > Security Options**
 - c. Double-click the following policy listed in the right-hand panel: **Network access: Sharing and security model for local accounts**
 - d. In the displayed dialog select the following item from the drop-down list: **Classic - local users authenticate as themselves**
- [MUST] 9. **Start > Settings > Updates and security > Windows Defender**:
- a. Click **Open Windows Defender Security Center**,
 - b. Click **App & browser control**, turn off **SmartScreen**:
 - SmartScreen for Microsoft Edge: **Off**
 - SmartScreen for Windows Store apps: **Off**
- [MUST] 10. **Date and Time** (Start > Windows System > Control Panel): Choose the time zone of your machine's location.
- [MUST] 11. **Network and Sharing Center** (Start > Windows System > Control Panel):
- a. Select **Change adapter settings**. Right-click the current active Ethernet, select **Properties > Configure**.
 - b. On the **Power Management** tab, clear all check boxes.



- [MUST] 12. **Programs and Features** (Start > Windows System > Control Panel):
- Click **Turn Windows features on or off**.
 - Enable .NET 3.5 by selecting the **.NET Framework 3.5 (includes .NET 2.0 and 3.0)** check box. This option requires an internet connection.
Note: If this procedure does not work as expected, or the computer has no internet access, install .NET 3.5 from the Windows installation media (see details for Windows 10 under <https://support.microsoft.com/en-us/kb/2734782>). If you do not have installation media, create them as described under <https://www.microsoft.com/en-us/software-download/windows10>.
 - To make sure that all the net.tcp components are properly initialized, Non-HTTP activation must be enabled. Expand the **.NET Framework 3.5 (includes .NET 2.0 and 3.0)** node and select the **Windows Communication Foundation Non-HTTP Activation** check box.
 - Select the **.NET Framework 4.7 Advanced Services** check box. Use the default values for sub items.
 - Select the **Internet Explorer 11** check box.
 - Select the **Telnet Client** check box.
 - Select the **TFTP Client** check box.
 - Reboot the PC if necessary.
- [MUST] 13. **Start > Settings > Apps > Default Apps:** For best user experience in OpenLAB Help and Learning, select Internet Explorer or Google Chrome as default **Web browser**.
- [MUST] 14. Disable **Compatibility View** in Internet Explorer.
- Start > Windows Accessories**, open Internet Explorer.
 - Click the Tools menu, and then click **Compatibility View Settings**.
 - Clear the **Display intranet sites in Compatibility View** check box.
- [MUST] 15. Enable the **Navigation pane**.
Start > Windows System, open **File Explorer**, then select **View > Navigation pane** from the ribbon and make sure that **Navigation pane** is selected.
- [MUST] 16. Disable **Windows Update service**
Start > Windows Administrative Tools > Component Services, select **Console Root > Services (Local)**, double click **Windows Update**, on **General** tab, set **Startup type: Disabled**, click **OK**.
- [PERFORMANCE] 17. **System** (Start > Windows System > Control Panel): Change performance options.
- Click **Advanced system settings**.
 - On the **Advanced** tab > **Performance** click **Settings**.
 - On the **Visual Effects** tab, select **Custom**.
 - Under **Custom**, select the following check boxes for better usability:
 - Smooth edges of screen fonts**
 - Show shadows under mouse pointer**
 - Show shadows under windows**
 - On the **Advanced** tab, select **Virtual Memory**: For optimum performance use the **Change** button to adjust the paging file size to a value of 2 to 3 times of the physical RAM on the PC. If possible, locate the paging file on a drive different from the system installation drive.
 - On the **Data Execution Prevention** tab, select **Turn on DEP for essential Windows programs and services only**.



[PERFORMANCE] 18. **System** (Start > Windows System > Control Panel): Change system properties.

- a. Click **Advanced system settings**.
- b. On the **Advanced** tab > **Startup and Recovery**, click **Settings**.
 - **System startup** section:
Change both **Time to display ...** fields from **30** to **3** sec.
 - **System failure** section:
Select **Automatically restart**, in the **Write debugging information** section
Select **Kernel memory dump** from the drop-down list.
- c. On the **System Protection** tab, in the **Protection Settings** section
Make sure that **Protection** for all drives is turned off. If required, click **Configure** and select **Disable system protection**.
- d. On the **Remote** tab, in the **Remote Assistance** section
Clear the check box **Allow Remote Assistance connections to this computer**.

[PERFORMANCE] 19. **Start > Settings > Personalization > Colors**: Turn **Transparency effects** off.

[OPTIONAL] 20. **Start > Settings > Personalization**:

- a. On the **Lock screen** page:
 - Under **Background**, select **Picture** or **Slideshow**.
 - Turn off **Get fun facts, tips, tricks, and more on your lock screen**.
 - Turn off **Show lock screen background picture on the sign-in screen**.
- b. On the **Start** page:
Turn off **Occasionally show suggestions in Start**.

[OPTIONAL] 21. **Start > Settings > Privacy**:

- a. On the **General > Change Privacy Options** page, turn off the following:
 - **Let apps use advertising ID to make ads more interesting to you based on your app usage (turning this off will reset your ID)**
 - **Let website provide locally relevant content by assessing my language list**
 - **Let Windows track app launches to improve Start and search results**
- b. On the **Location** page, make sure **Location** for this device is off. If not, click **Change** to turn it off.

[OPTIONAL] 22. Right click on **Start**, select **Run** from the context menu, type **gpedit.msc** in the **Run** field:

- a. Navigate to **Local Computer Policy > Computer Configuration > Administrative Templates > System > Logon**.
- b. Set **Don't display the Getting Started welcome screen at logon** to **Enabled**.

[OPTIONAL] 23. **Recycle Bin Properties**: (right-click on desktop icon **Recycle Bin**) Select the following options:

- a. **Custom size**: Select a size corresponding to approximately 10% of the complete disk space for the drive.
- b. Select **Display delete confirmation dialog**.

Repeat these steps for all drives of your computer.

[OPTIONAL] 24. **Region** (Start > Windows System > Control Panel): Language for non-Unicode programs:

On the **Administrative** tab, click **Change system locale...** From the dropdown list, select **English (United States)**.

Note: Do not change system locale if you are using an **English, Portuguese, Japanese** or **Chinese** Operating System.



[OPTIONAL] 25. **Start > Settings > Personalization.** On the **Taskbar** page, under **Combine taskbar buttons** select **When taskbar is full.**

This will simplify switching between open CDS instances.