

AutoCONFIG scroll bar issue created by a Windows update on computers with version 1909

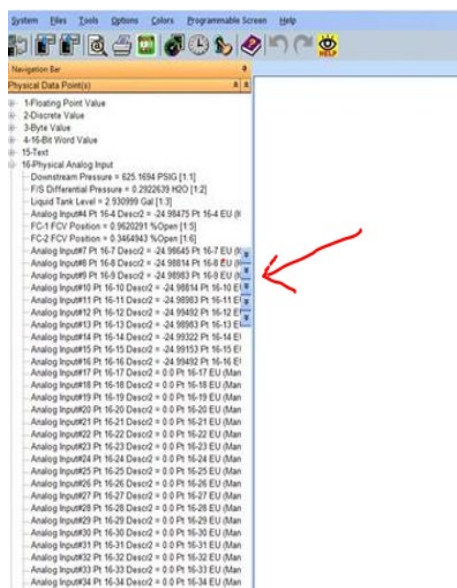
Key Words

AutoCONFIG, Windows Update, Scroll bar

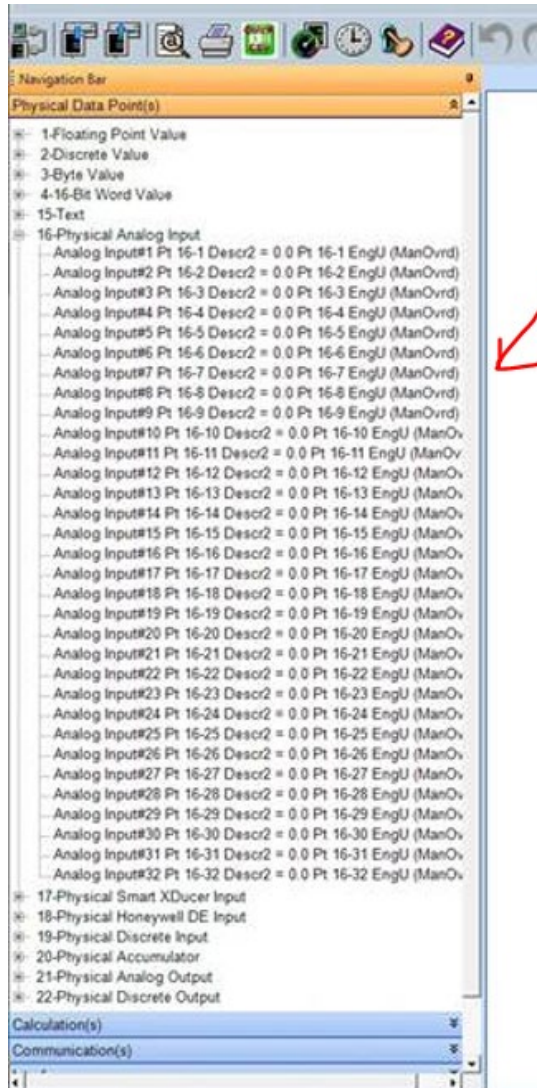
Introduction

The problem has been identified that the scroll bars within the navigation bars on the left side of the AutoCONFIG window have been disappearing without any changes on users' program. Upon investigation, it has been found that the issue is a consequence of a Windows update in the device with the AutoCONFIG program. This tech bulletin provides instructions on how this issue can be resolved.

If you are encountering the issue, the AutoCONFIG window will look as shown below.



However, under normal conditions the same screen should look as shown below with the scroll bar.

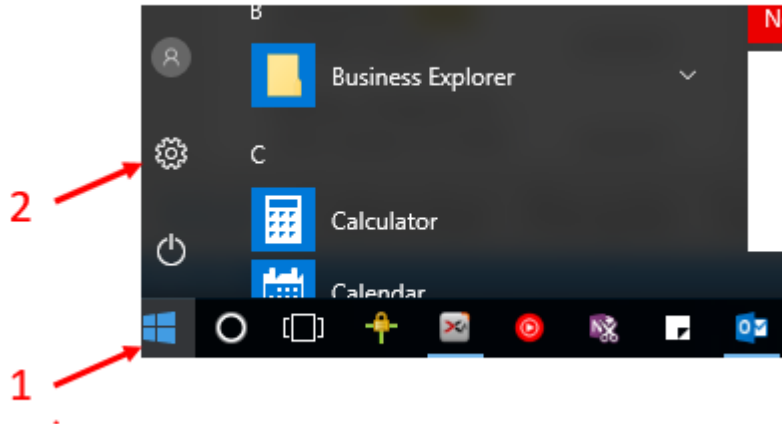


Solution

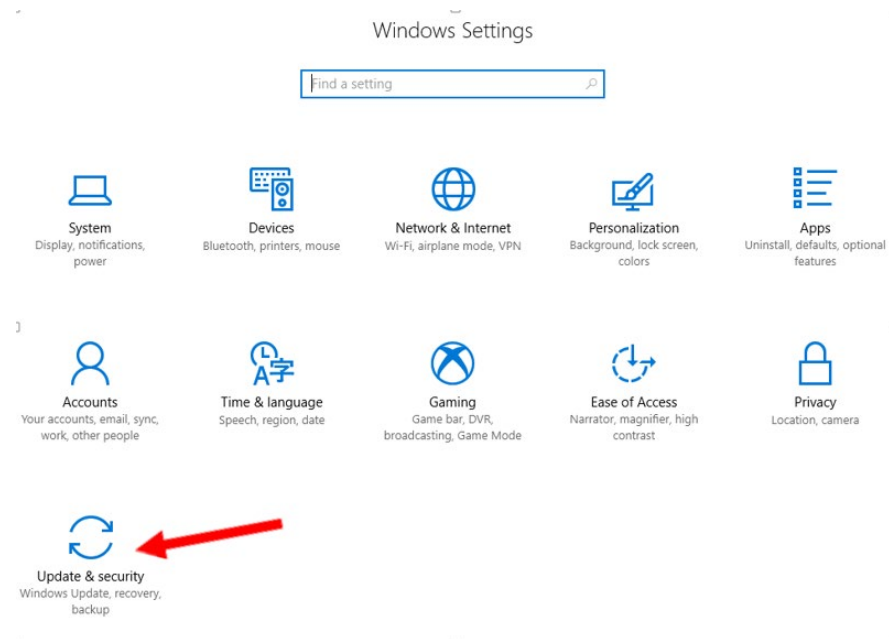
This issue can be resolved simply by updating the windows device to the most recent updates. To get the next version of the update that would replace the problematic version with KB5001337 security update, simply go to the settings in your device, then click on Update and Security.

When you are on that page, click “Check for Updates” and download the latest updates that are suggested for your device.

Step 1.



Step 2.



Step 3.

Windows Update

Update status



Your device is up to date. Last checked: Today, 12:45 PM

Check for updates



[Update history](#)

USA

27 Forge Parkway

Franklin, MA 02038

Ph: (713) 272-0404

orders.process.us@thermofisher.com

Find out more at thermofisher.com/autoconfig

© 2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.