

Prometheus Control PC

This document specifies the system configuration required for Prometheus systems (NT.48, NT.Plex, and Panta), as well as the configuration of PCs currently delivered with such instruments.

Systems connected to an NT.Robotic Autosampler have additional requirements listed in a separate document.

System requirements for Prometheus control PCs

Device Type	Prometheus NT.48 or NT.Plex	Prometheus Panta
Software	PR.ThermControl PR.ChemControl PR.TimeControl PR.Stability Analysis PR.ThermControl-CFR	PR.Panta Control PR.Panta Analysis
.NET Framework	.NET Framework 4.5.1 or higher .NET Framework 4.5.1 is included in the installer	.NET Framework 4.6.2 or higher .NET Framework 4.6.2 is included in the installer
Windows version	Windows 10 (x64) with Anniversary Update Windows 11	
Windows updates	Should not be installed during measurements	
Virus scanner	Tested with Windows Defender	
Ports	1x 100 Mbit RJ45 Ethernet Port (Device) 2x USB for Ethernet adapter (Mouse, Network adapter)	
Network config	Internal Ethernet Port must use IP address 192.168.1.1 and subnet mask 255.255.255.0.	Internal Ethernet Port must use IP address 172.16.4.3 and subnet mask 255.255.255.0.
	Connection to corporate network via USB to RJ45. PC can be added to domain if domain settings (firewall, updates) are supported.	
Firewall settings	Avoid firewalls between control PC and device. If a firewall is present, please allow the following connections:	
	Port 12000/9020, TCP/UDP, In/Out; Port 137/3702/5355, UDP, In/Out; ICMPv4/ICMPv6, In/Out	Port 5001, TCP, In/Out; Port 8888, TCP, In/Out; Port 22, TCP/UDP, In/Out; Port 9090, TCP, In/Out; Port 2318, TCP/UDP, In/Out
Display	Full HD @ 125%	
Memory	8 GB or higher	

CPU	Intel Core i5, 6th generation or higher	Intel Core i5, 8th generation or higher
Hard drive	20 GB available space, SSD recommended	
Administrator privileges	Only required for installation	

Important notes

- The Ethernet port which is directly connected to the Prometheus instrument (it must be a direct port-to-port connection, do not use a connection via switch or hub) must have the static IP-Address **192.168.1.1** for Prometheus NT.48 and Prometheus NT.Plex, or **172.16.4.3** for Prometheus Panta, and the subnet mask **255.255.255.0**. Additionally, the port **9020** must not be blocked.
- Avoid firewalls, they can harm the real-time data transfer between the Prometheus instrument and the control PC.
- The software was tested to function with Windows Defender. Other virus scanners might harm the real-time data transfer between the Prometheus instrument and the control PC.
- Please install TeamViewer (TeamViewer GmbH) to allow NanoTemper Technologies to support you via remote control.
- An external computer mouse is necessary to access all software features.