

Monolith NT.115 Control PC Requirements

MO.Control is the latest software from NanoTemper Technologies for controlling your Monolith NT.115, Monolith NT.115Pico, Monolith NT.LabelFree and Monolith NT.Automated system. MO.Control has replaced the old software NT.Control for most Monolith systems*.

System configuration needed for NT.115 control PCs

Device Type	NT.115, NT.115 Pico, NT.LabelFree
Installed software	MO.Control MO.Affinity Analysis
.NET Framework	.NET Framework 4.5.1 or higher, .NET Framework 4.5.1 is included in the installer
Windows version	Windows 7/10/11 (x86 or x64)
Windows updates	Should not be installed during measurements
Virus scanner	Tested with Windows Defender
Ports	1x 100 Mbit RJ45 Ethernet Port (Device), 3x USB (Mouse, data stick, network)
Network config	Internal Ethernet Port must use IP address 192.168.1.1 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
Additional requirements	%ProgramFiles(x86)\CommonFiles\NanoTemper\Control\Config.xml needs to be copied from NTT supplied computer to own computer
Display	Full HD @ 125%
Memory	8 GB or higher
CPU	Intel Core i5, 6th 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 Monolith 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** and the Subnetmask **255.255.255.0**. Additionally, the port **9020** must not be blocked.
- Avoid firewalls, they can harm the real-time data transfer between the Monolith 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 Monolith instrument and the control PC.
- Please install TeamViewer (TeamViewer GmbH) to allow NanoTemper Technologies to support you via remote control.
- %ProgramFiles(x86)\CommonFiles\NanoTemper\Control\Config.xml must be copied to the operating computer