

Modern instrument, familiar computer

Details on the Nicolet Summit integrated computer



Unlike other laboratory equipment that uses proprietary integrated computers, the Thermo Scientific™ Nicolet™ Summit FTIR Spectrometer includes a standard Windows® 10 computer that acts just like a common laptop. The instrument can easily be added to the company domain and equipped with company security policies like:

- Access control policies (ACPs)
- Change management policies (CMPs)
- Information security policies (ISPs)

These system protections will be automatically deployed on the Nicolet Summit Spectrometer once it is on the domain, making it ideal for use in regulated environments.

Advantages of an integrated computer:

- 1 Save money** – no separate PC required
- 2 Full Wi-Fi connectivity** – connect to networks and send data directly to other devices
- 3 Seamlessly switch interfaces** – go from touchscreen to desktop software interface
- 4 Powerful capabilities** – automated background collection (Smart Background); LightBar visual results
- 5 Save bench space** – smaller footprint for more samples and better portability

Features	Summit	Summit Pro
Integrated Windows 10 computer with 64 GB eMMC storage	✓	✓
Free access to Thermo Fisher Cloud with 1 TB storage	✓	✓
9th Generation Intel® GPU	✓	✓
DRAM	4 GB	8 GB
Processor	Intel® Celeron® Dual Core	Intel® Pentium® Quad Core

Flexible connectivity and data storage

The data collected on the Nicolet Summit Spectrometer is stored in a database. The default database, MariaDB®, resides on the instrument itself, but a variety of other databases can also be used. With the Remote Database function, you can also configure the instrument to store data in another database location, like a secured company server (local or networked database locations).

Designed for long-life

The Nicolet Summit Windows 10 computer has been specifically designed for extended use in a laboratory environment. It includes:

1. eMMC hard-drive with no moving parts
2. Conformal electronics coating to protect against chemicals, dust, and temperature swings
3. Next-generation Intel processor, supported for >15 years
4. Components from Intel's 'Internet of Things' (IoT) long-term product roadmap
5. Common form-factor computer module to keep replacement costs low

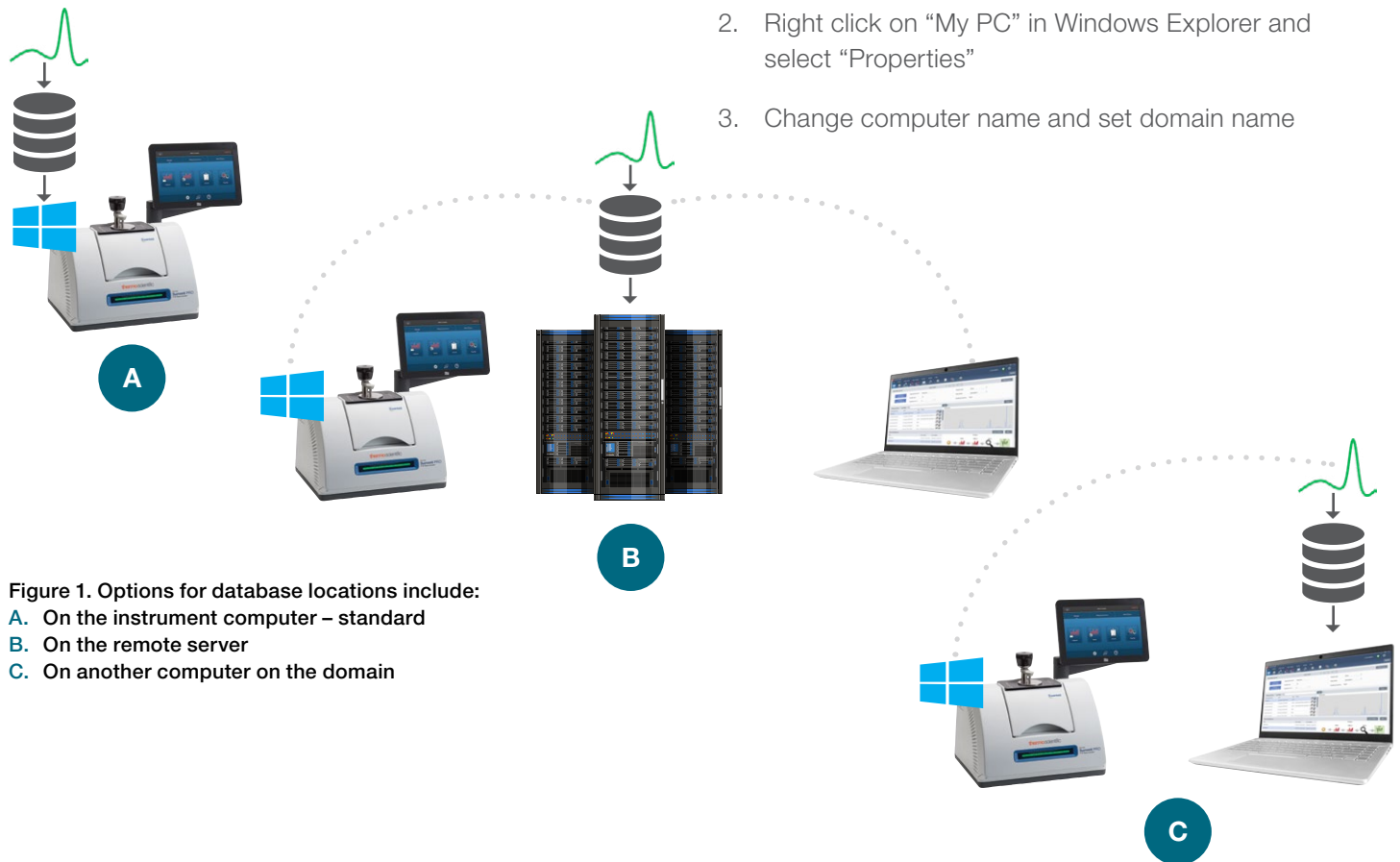


Figure 1. Options for database locations include:

- A. On the instrument computer – standard
- B. On the remote server
- C. On another computer on the domain

Database	Supported versions
AuroraDB	MySQL 5.6
	MySQL 5.7
MariaDB	10.0
	10.1
	10.2
	10.3
Oracle	11g Release 2
	12c Release 1
	12c Release 2
	18c
SQLServer	2012
	2014
	2016
	2017

How to connect the spectrometer to a company domain

1. An IT administrator logs into computer with domain-add permission
2. Right click on "My PC" in Windows Explorer and select "Properties"
3. Change computer name and set domain name

Find out more at thermofisher.com/summit