



# Agilent OpenLAB ECM Software and Hardware Requirements

For Revision 3.5

The table below lists the software and hardware requirements for Agilent OpenLAB ECM 3.5. This applies to OpenLAB ECM Enterprise and OpenLAB ECM Workgroup configurations.

## Software and Operating System Requirements

		Compatible Version	
Software Application Modules	Related Components	OpenLAB ECM Server	Version 3.5
		OpenLAB ECM Options CD	Version 3.5
System Requirements	ECM Server	Operating System*	Windows 2012 Server R2 (64 bit) US/English version
		Regional Settings	English US
		.Net Framework	Version 3.5, Version 4.5.1
		Browser	Internet Explorer 11.x (32 bit IE only)
	ECM Client	Operating System*	Windows 7 (32 or 64 bit) Windows 10 (32 or 64 bit) Windows 2012 Server R2 (64 bit)
		Regional Settings	English US or Western European. English US is required for client machines running Schedulers or any add-ons that generate XML data or perform filtering.
		.Net Framework	Version 3.5, Version 4.5.1
		Adobe Acrobat	Version 10, Version 11
		MS Office	Version 2010, Version 2013 (32 bit Office only)
		Browser	Internet Explorer 11. x (32 bit IE only)
	Database Server	Oracle Version	Oracle 12c Server (64 bit) and Oracle 11g R2 Client (32 bit, version 11.2.0.4) Standard or Enterprise Edition
		SQL Server Version**	SQL Server 2012 SP2 or SP3 Standard or Enterprise SQL Server 2014 Standard or Enterprise (Collation setting of SQL_Latin1_General_CP1_CI_AS is required)
		Regional Settings**	English US.

\* The US/English version of the Operating System is required for ECM Servers and ECM Schedulers

\*\* If the OS settings DURING INSTALLATION OF THE OPERATING SYSTEM are for US English then the default installation settings for SQL Server will be appropriate. Registry value changes will otherwise be necessary. See the ECM Installation Guide for details. Running a database server on a virtual machine is not a recommended configuration.



**Agilent Technologies**

## Hardware Requirements

### OpenLAB ECM Web Server Minimum Hardware Requirements\*

Component	Description	Notes
Processor	2.0 Ghz or higher	Dual Core or dual processor recommended
Memory	4 GB	8 GB recommended. Varies based upon concurrent users.
Hard Drive	100 GB	
Optical Drive	CDROM Drive	Recommended for software installation
Network Interface	100 Mbps	

\* Or comparable Virtual Machine using VMWare ESX 5.x or higher.

### OpenLAB ECM File Transfer Server Minimum Hardware Requirements

Component	Description	Notes
Processor	2.0 Ghz or higher	Dual Core or dual processor recommended
Memory	4 GB	8 GB recommended
Hard Drive	500 GB	A dedicated 100 GB partition is needed for the FT Cache
Optical Drive	CDROM Drive	Recommended for software installation
Network Interface	100 Mbps	

### OpenLAB ECM Application Server Minimum Hardware Requirements

Component	Description	Notes
Processor	2.0 Ghz or higher	Dual Core or dual processor recommended
Memory	4 GB	8 GB or more recommended
Hard Drive	500 GB	At least 50 GB of free disk space is needed for temporary file processing activities
Optical Drive	CDROM Drive	Recommended for software installation
Network Interface	100 Mbps	

### SQL Database Server Minimum Hardware Requirements

Component	Description	Notes
Processor	2.0 Ghz or higher	Dual Core or dual processor recommended
Memory	8 GB	16 GB recommended
Hard Drive	100 GB	Multiple drives are recommended for added performance
Optical Drive	CDROM Drive	Recommended for software installation
Network Interface	100 Mbps	

## OpenLAB ECM Client Minimum Hardware Requirements

Component	Description	Notes
Processor	1.5 GHz or higher	Dual Core or dual processor recommended
Memory	4 GB for Windows 7 8 GB for Windows 10 8 GB for Windows Server 2012 R2	
Hard Drive	250 GB	
Optical Drive	CDROM Drive	Recommended for software installation
Network Interface	10 Mbps	100 Mbps recommended
Graphics	1024x768 minimum resolution	1280x1024 or higher resolution recommended

\* Or comparable Virtual Machine using VMWare ESX 5.x or higher.

## OpenLAB ECM Scheduler System Minimum Hardware Requirements

Component	Description	Notes
Processor	2.5 GHz or higher	Dual Core or Quad Core recommended
Memory	4 GB for Windows 7 8 GB for Windows 10 8 GB for Windows Server 2012 R2	
Hard Drive	500 GB	Drive used for the scheduler inbox should be larger than amount of files being uploaded.  The Scheduler needs disk space available on the drive designated by the Temp variable. To process SSZip files, at least 3x the maximum size of an SSZip archive is needed.
Optical Drive	CDROM Drive	Recommended for software installation
Network Interface	100 Mbps	
Graphics	1024x768 minimum resolution	1280x1024 or higher resolution recommended

It is not recommended to run the Scheduler on the same machine that is hosting the OpenLAB ECM Server.

## All-in-One OpenLAB ECM Server or Workgroup Server Minimum Hardware Requirements

Component	Description	Notes
Processor	4.0 Ghz or higher	Dual Core or dual processor recommended
Memory	12 GB	24 GB or higher recommended. If the database is also running on the same machine, review the recommendations from Oracle or Microsoft as well.
Hard Drive	2 TB	Disk space sufficient for 3 years of file storage is recommended.
Optical Drive	CDROM Drive	Recommended for software installation
Network Interface	100 Mbps	1 GB Mbps recommended

[www.agilent.com/chem/openlab](http://www.agilent.com/chem/openlab)

© Agilent Technologies, Inc., 2016  
Published in USA, August 30, 2016  
Publication Number 5991-6503EN



**Agilent Technologies**