

# Installing Polara RS

This document describes how to install Polara RS on a PC. Full details about installing on a lab system computer and an offline computer are in the *Polara RS Lab System User Guide* that is available after you complete this installation.

This document contains the following information:

- “Installation Requirements” on page 2.
- “Contents of the Installation Key” on page 3.
- “Installing on a Computer” on page 4.

## **Installation Requirements**

The lab computer requires the following hardware and software to run Polara RS.

- Computer Software
  - Microsoft Windows XP with the latest service packs and updates installed
  - Microsoft Internet Explorer 5.01 or later
  - Microsoft .NET Framework 1.1 Service Pack 1 and Framework 2.0 (supplied)
- Computer Hardware
  - 90 Mhz Intel Pentium-class processor (minimum requirement, at least 200 Mhz recommended)
  - 32 MB of RAM or more (at least 96 MB recommended)
  - 150 MB of hard disk storage for software, plus storage space for run-time data (at least 200 MB recommended)
  - 1 USB port
  - COM ports - one for the RapidStak and one for each instrument
  - 800 x 600 pixel or higher resolution video display

## Contents of the Installation Key

The Installation Key, which is a storage device, has files organized in directories according to their purpose.

Files and Directories on the Installation Key		Contents	Description
Files	ReadMe.htm		Links to this document and Release Notes.
	Installing Polara RS.htm		This document and its style sheet.
	Installing Polara RS.css		
	Release Notes.txt		Release notes for Polara RS.
Directories	Setup	CmRuntimeUser.exe PolaraRSSetup.msi	Files for installing Polara RS and enabling it for lab system use.
	DotNet	dotnetfx1.1.exe dotnetfx2.0.exe	Microsoft installation files required for Polara RS.
	Firmware	bootManager.exe firmware.hex	RapidStak files for updating firmware.
	Drivers	*.msi	Installation files for each driver except the RapidStak.
	Adobe	AdbeRdr707_en_US.exe	Installation file for Acrobat Reader

Read the Release notes to learn about Polara RS and driver issues that are not documented elsewhere.

## Installing on a Computer

To install Polara RS on a lab computer, ensure you meet all the requirements in “[Installation Requirements](#)” on [page 2](#). When you are ready to install, follow the procedure “[To install Polara RS](#)” on [page 4](#).

**Note** To run assays on the lab system computer with instruments in normal mode, you must plug the Installation Key into a USB port.

### To install Polara RS

1. Use Windows Explorer to access the root directory on the Installation Key.
2. Install Microsoft’s .NET Framework, as follows:
  - a. Open the DotNet directory.
  - b. Double click **dotnetfx1.1.exe** and complete the installation procedure.
  - c. Double click **dotnetfx2.0.exe** and complete the installation procedure.
  - d. Return to the root directory.
3. Install Adobe’s Acrobat Reader, if it is not installed, as follows:
  - a. Open the Adobe directory.
  - b. Double click **AdbeRdr707\_en\_US.exe** to start the installation wizard. Complete the wizard.
  - c. Return to the root directory.
4. Install Polara RS as follows:
  - a. Open the Setup directory.
  - b. Double click **CMRuntimeUser.exe** to start the CodeMeter installation wizard. Complete the wizard.
  - c. Double click **PolaraRSSetup.exe** to start the Polara RS installation wizard. Complete the wizard.
  - d. Return to the root directory.
5. Install drivers as follows:

- a. Open the Drivers directory.
  - b. For each instrument, double click its **.msi** file to start the driver installation wizard for the instrument. Complete each wizard.
6. If the computer is not a lab system computer, remove the Installation Key and plug it into the lab system computer.
  7. If the computer is a lab system computer, contact your IT department to ensure your computer is not using Hyper-Threading. When Hyper-Threading is enabled on a computer, Polara RS runs correctly, but execution can slow significantly.