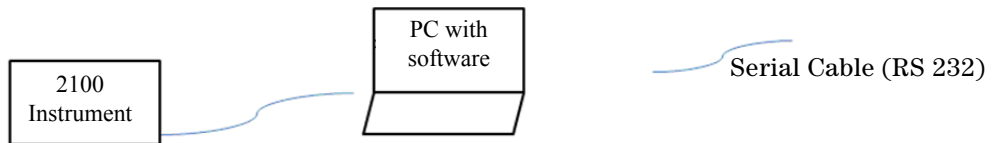


## Agilent 2100 Bioanalyzer – PC connection instructions

This document provides information about the Agilent 2100 Bioanalyzer and its options for connection to PC's. There are two options depending on what type of communication ports are available for the existing PC.

### Instrument connection to PC's via serial (COM) ports Recommended, standard connection - all drivers are preinstalled

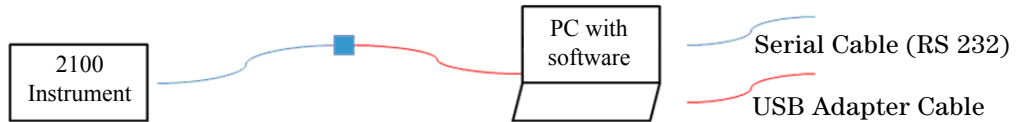


- 1 Connect 2100 Bioanalyzer instrument via Serial cable to PC
- 2 Turn on 2100 Bioanalyzer instrument and PC
- 3 Start 2100 Expert Software and choose correct COM port



## Instrument connection to PC's via USB ports In case no serial port is available

**Agilent bundle Laptop - all drivers are preinstalled  
(Concerns only product G2953CA)**



- 1 Connect 2100 Bioanalyzer instrument via serial cable and adaptor cable to PC
- 2 Turn on 2100 Bioanalyzer instrument and PC
- 3 Start 2100 Expert Software and choose correct COM port

### External PC's

**Requires driver installation from 2100 Expert Software CD-ROM !**



- 1 Install USB-Adapter driver from 2100 Expert Software CD-ROM, or download the driver online
  - a Insert Agilent CD-ROM, then choose directory and file :  
  \Support\Driver\5188-8031\CDM20824\_Setup.exe
  - b Start file CDM20824\_Setup.exe
  - c Reboot PC
- 2 Connect 2100 Bioanalyzer instrument via Serial cable and Adaptor cable to PC
- 3 Turn on 2100 Bioanalyzer instrument
- 4 Start 2100 Expert Software and choose correct COM port



G2938-90125 Rev. B  
Edition: 02/2017

For Research Use Only  
Not for use in diagnostic procedures

Printed in Germany

© Agilent Technologies 2017

Agilent Technologies  
Hewlett-Packard-Strasse 8  
76337 Waldbronn, Germany