

October 11, 2013

---

## Table of Contents

---

1. What's New?
    - 1.1 2100 Bioanalyzer Support Tool Box, version 1.
  2. System Configuration
    - 2.1 System Requirements
  3. Software Installation and Setup
    - 3.1 Installation Instructions
    - 3.2 How to Use
    - 3.3 Remove Program Instructions
  4. Known Issues
    - 4.1 Buttons will not launch PDF files
  5. Navigation Guide
- 

### 1. What's New?

---

#### 1.1 2100 Bioanalyzer Support Tool Box, version 1.

The 2100 Bioanalyzer Support Tool Box (SuTB) provides content to maximize the efficiency of the Agilent 2100 Bioanalyzer System. Content includes how-to familiarization videos, maintenance guidelines, essential measurement practices, assay conditions, solutions to common issues, FAQs, and manuals. The SuTB contains primarily off-line content to allow access to documents without the need for an internet connection.

---

### 2. System Configuration

---

#### 2.1 System Requirements

The 2100 Bioanalyzer SuTB was tested on HP PCs with Windows 7. Other requirements specified for viewing are:

##### 2.1.1 Flash

Flash Player 10 or later (<http://www.adobe.com/go/getflash>), and one of the following browsers:

- Windows: Internet Explorer 6 and later, Firefox 1.x and later, Safari 3 and later, Google Chrome, Opera 9.5 and later

##### 2.1.2 PDF viewer

- Agilent supports Adobe Acrobat Reader X. It may be downloaded directly from Adobe (<http://get.adobe.com/reader>).
- 

### 3. Software Use

---

#### 3.1 Installation Instructions

- Double-click **Agilent 2100 Bioanalyzer Support Tool Box\_Setup.exe**. Follow prompts for installation.

*NOTE: Windows 7 users must have administrator rights.*

#### 3.2 How to Use

- To open the program, click on: Start menu > All Programs > Agilent 2100 Bioanalyzer Support Tool Box folder > Agilent 2100 Bioanalyzer Support Tool Box. Click on buttons within the program to access information.

### 3.3 Remove Program Instructions

- Click on: Start menu > All Programs > Agilent 2100 Bioanalyzer Support Tool Box folder > Uninstall. Follow prompts to remove the program.

## 4. Known Issues

### 4.1 Buttons will not launch PDF files

This is often caused by security settings in Adobe. Open Adobe Acrobat Reader X > Edit > Preferences > General > Uncheck "Enable Protected Mode as startup".

## 5. Navigation Guide

