

# How to Download, Install, and Register RTCA Software Pro 2.8.1

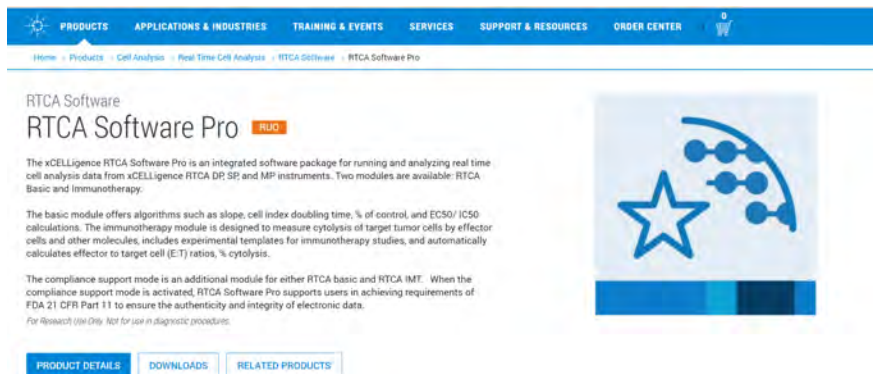
## Compatible instruments:

This software is compatible with the following xCELLigence Real-Time Cell Analysis (RTCA) models:

- xCELLigence RTCA SP
- xCELLigence RTCA DP
- xCELLigence RTCA MP
- xCELLigence RTCA TP\*
- xCELLigence RTCA HT
- xCELLigence RTCA S16
- xCELLigence RTCA iCELLigence\*\*

The latest version of the software can be downloaded from our website:

<https://www.agilent.com/en/product/cell-analysis/real-time-cell-analysis/rca-software/rca-software-pro-741236>



PRODUCTS APPLICATIONS & INDUSTRIES TRAINING & EVENTS SERVICES SUPPORT & RESOURCES ORDER CENTER

Home > Products > Cell Analysis > Real Time Cell Analysis > RTCA Software > RTCA Software Pro

### RTCA Software

## RTCA Software Pro BUY

The xCELLigence RTCA Software Pro is an integrated software package for running and analyzing real time cell analysis data from xCELLigence RTCA DP, SP, and MP instruments. Two modules are available: RTCA Basic and Immunotherapy.

The basic module offers algorithms such as slope, cell index doubling time, % of control, and EC50/IC50 calculations. The immunotherapy module is designed to measure cytotoxicity of target tumor cells by effector cells and other molecules, includes experimental templates for immunotherapy studies, and automatically calculates effector to target cell (E:T) ratios, % cytotoxicity.

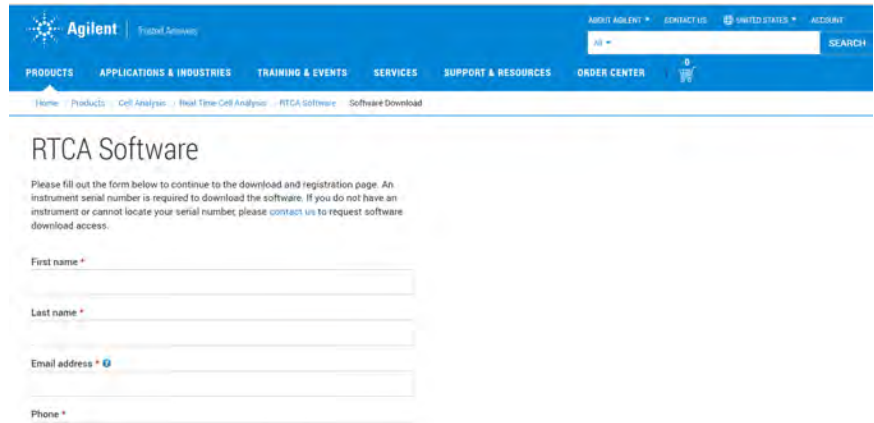
The compliance support mode is an additional module for either RTCA basic and RTCA IMT. When the compliance support mode is activated, RTCA Software Pro supports users in achieving requirements of FDA 21 CFR Part 11 to ensure the authenticity and integrity of electronic data.

For Research Use Only. Not for use in diagnostic procedures.

PRODUCT DETAILS DOWNLOADS RELATED PRODUCTS

1. Select "Downloads"

You will be routed to the page shown on the right. If you have previously downloaded software, skip to step number four.



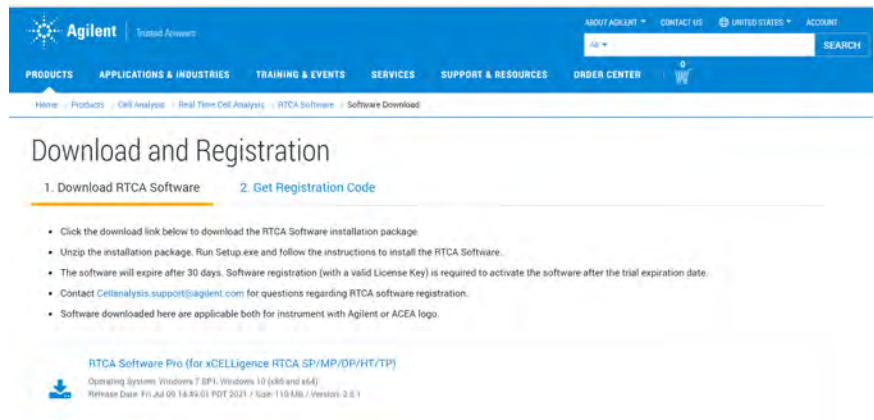
2. Fill out the form. When prompted for the serial number (SN), enter the SN listed on the instrument that operates inside the incubator (with cradles for plates).

The SN can be located on the back of the instrument. Use the list below to ensure you have the correct serial number:

- xCELLigence RTCA SP: Single plate (1 x 96-well plate); SN will start with 26-.
- xCELLigence RTCA MP: Multiple plates (6 x 96-well plates); SN will start with 29-.
- xCELLigence RTCA DP: Dual plate (3 x 16-well plates); SN will start with 32-.
- xCELLigence RTCA HT: High-throughput model (384-well plate); SN will start with 34-.
- xCELLigence RTCA S16: Pilot scale model (1 x 16-well plate); SN will start with 58-.
- xCELLigence RTCA TP\*: Only supported in parts of Asia.
- xCELLigence RTCA iCELLigence\*\*: (3 x 8-well plates); No longer supported by Agilent.

3. Select "Next" at the bottom of the page.

You will be directed to download the software. Please make sure that the name of the software on the next page is the correct one you intend to download.



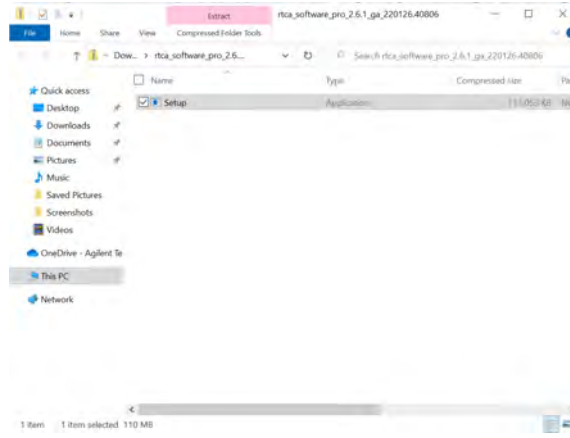
4. Click on the software link and follow the wizard.

- Once you download the software from the Agilent website, you will be unable to download a second software unless you delete your cookies or open a window in "private" or "incognito" mode in your browser. Select "New InPrivate window" in Microsoft Edge or "New Incognito window" in Google Chrome.
- Copy and paste the following link into a private window: <https://www.agilent.com/en/product/cell-analysis/real-time-cell-analysis/rtca-software/rtca-software-pro-741236>
- Return to step 2.

The download will appear as a pop-up under "Downloads" and will be available in the "Downloads" folder.



5. Once the software download is completed, select "Open file".



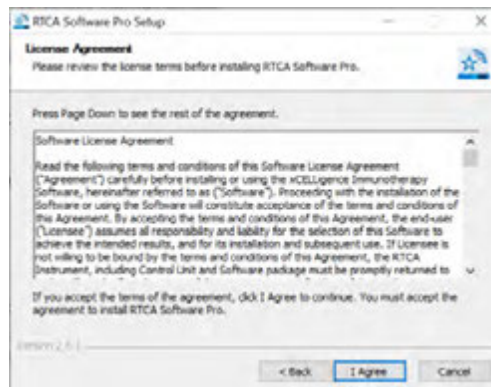
6. Place the zip file onto your desktop and extract all files.

7. Double click on the file titled "Setup.exe".

The RTCA Software Pro Setup Wizard window will open.

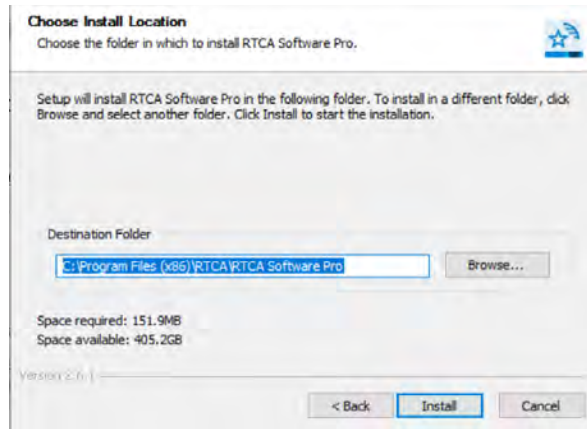


8. Click "Next".

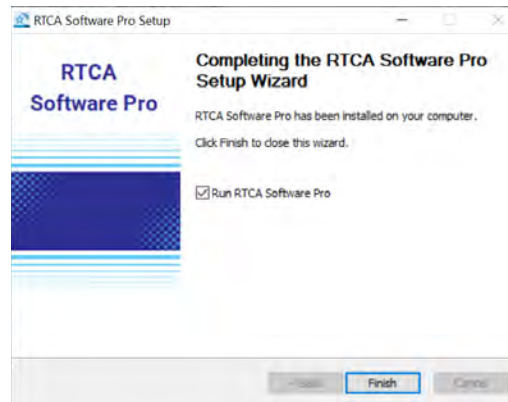


9. Review the software license agreement. If you accept the terms, click "I Agree".

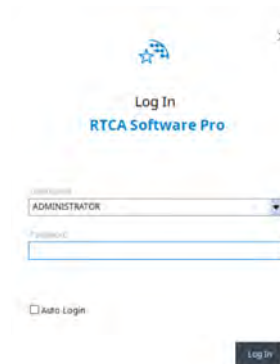
The RTCA Software Pro Setup Wizard  
(continued).



10. Choose which folder to install the software to. Click "Install".

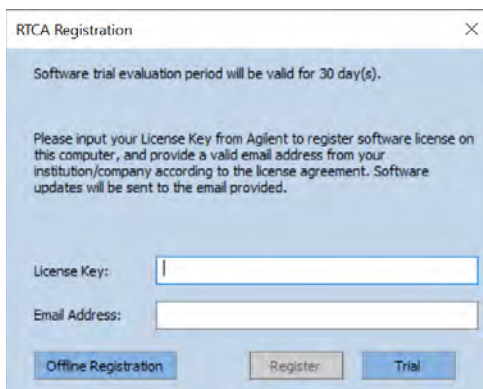


11. Click "Finish".



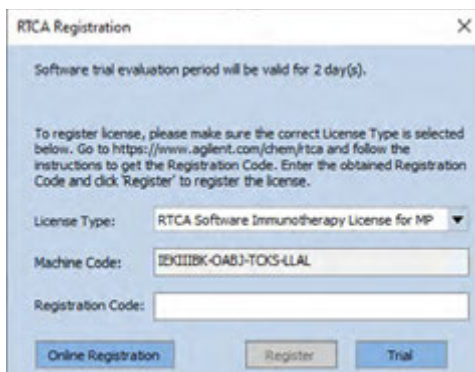
12. Log in to the software. The password is "administrator".

A pop-up window will appear before opening the software.



13. Enter your license key and email and click "Register". If you do not have a license key, select "Trial".

If the computer is not connected to the internet:

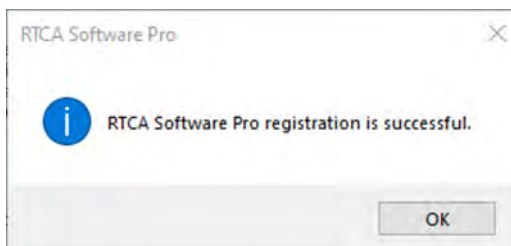


14. Click "Offline Registration" to switch to offline registration mode. Select the "License Type" you purchased from the drop-down menu and copy the "Machine Code" displayed in the window.

15. On a computer that is connected to the Internet, go to <https://www.agilent.com/en/product/cell-analysis/real-time-cell-analysis/rtca-software/rtca-software-pro-741236>:

- Fill out the form with the xCELLigence station serial number, then click "Next".
- On the "Download and Registration" page, click "Get Registration Code".
- Enter valid "Machine Code" and "License Key".
- Click "Get Registration Code".
- Copy the "Registration Code", enter it to the "Software Registration" window, and click "Register"

A message box will pop up, indicating the result of the registration.



NOTE: A valid email address of your company/institution must be provided according to the license agreement. Software update information will be sent to the email provided.

NOTE: A license key is required for software registration or license transfer. Keep a safe record of the license key assigned to each computer or user. You can continue to register other licenses after registering one license successfully.

[www.agilent.com/lifesciences/rtca](http://www.agilent.com/lifesciences/rtca)

DE40155023

This information is subject to change without notice.

© Agilent Technologies, Inc. 2024  
Published in the USA, June 18, 2024  
5994-7271EN

