

xPONENT 3.1 Upgrade Questions and Answers (Q&A)

Thank you for your interest in our Luminex **xPONENT 3.1** Software Solutions.

This Upgrade Q&A document will help ensure that your current Luminex System and PC meet the minimum specifications required to allow an upgrade to the new software.

The System Requirements Checklist will assist in determining that your system components are upgrade ready.

The Upgrade Order Checklist will ensure that your customized **xPONENT 3.1** Upgrade Kit contains your required items.

Please contact Luminex Technical Support if you need assistance in locating any of the requested information.

Luminex Technical Support is available to users in the U.S. and Canada by calling 1-877-785-BEAD (-2323). Users outside of the U.S. and Canada can contact us at +1-512-381-4397. Inquiries may also be sent by email to support@luminexcorp.com.

Using this information, we will work with your sales representative to ensure you have all the items required for a successful upgrade to our new software solution, **xPONENT 3.1**.

We look forward to hearing from you!

Thank you,
Luminex Corporation

System Requirements Checklist

Please record the following information. If you need assistance locating any particular item or if you have any questions, Luminex Technical Support will be happy to assist you.

1) Contact and location information:

Name: _____

Phone number: _____

Address 1: _____

Address 2: _____

City, State,
Zip: _____

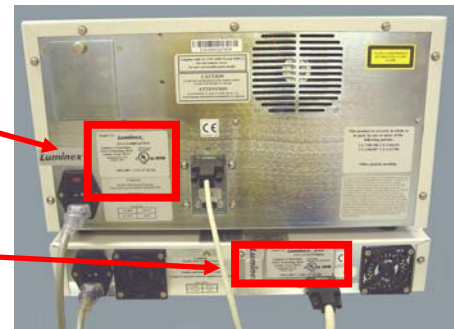
Partner through which you purchase Luminex systems: _____

2) Please provide the serial number of your Luminex System: LX100

This is located on the back of your LX100 or LX200. For both, the serial number begins with "LX100".

3) Please provide the serial number of your XY platform: LXY

This is located on the back of your XY platform. The serial number begins with "LXY".



4) If applicable, please provide your xPONENT 3.0 license code _____
This is located in the xPONENT 3.0 Admin Page → Licensing Tab

5) Please determine if your PC and monitor meet the following minimum specifications:

Operating System: Microsoft® Windows® XP Professional, SP3

Processor: 1.86 GHz Intel® Core™2 or higher (2.66 GHz recommended)

Memory: 2 G RAM or higher (4GB recommended)

Hard drive: 80 GB

External Drive: DVD

Screen resolution and number of colors: 1280 x 1024, 32 bits

To locate Operating System, processor and memory (RAM) specifications, click Start → (right-click) 'My Computer' → Properties. The information is listed on the 'General' tab that automatically appears.

To locate hard drive size and available hard drive space, click Start → My Computer → (right-click) Local Drive (C:) → Properties. 'Capacity' is size of the hard drive, and 'Free Space' is the amount of available hard drive space.

To locate resolution and number of colors, right click on an empty section of the desktop → Properties → Settings tab. If necessary, move the slider under 'Screen Resolution' to the right to determine if 1280 x 1024 is an option. The drop-down menu for 'Color Quality' displays the available numbers of colors.

6) Does your PC meet the requirements listed above? Yes No

7) Does your monitor have the capability to display 1280 x 1024? Yes No

Upgrade Order Checklist

Please provide the following information for your customized **xPONENT 3.1** Upgrade Kit.

- 1) Is the LX200 with xPONENT 3.1 being used or will be used in a clinical/ diagnostic lab? Yes No
This will help determine the modules you will or will not need.

- 2) From what version of Luminex software are you upgrading?
 LDS 1.7 SP1 LDS 1.7 SP2 IS 2.1 IS 2.2 IS 2.3 xPONENT Other: _____
To determine which Luminex software version you are using, launch the Luminex software. The information is displayed on the splash screen. It can be also displayed once the Luminex software is running by clicking Help → About.

- 3) Would you like to order a new PC? Yes No
*Please note that if your PC does not meet the minimum specifications, you will need to purchase a PC meeting those specifications in order to use **xPONENT 3.1** software.*

- 4) Would you like to order a new monitor? Yes No
If so, please select one of the following options:
 Color touch screen monitor
 Color flat panel monitor
*Please note that if your monitor does not meet the minimum specifications, you will need to purchase a monitor with those specifications in order to use **xPONENT 3.1** software.*

- 5) Do you require the Basic Security Module? Yes No
The security module provides security through user login and passwords. It is automatically included when you purchase the 21CFR Part 11 module.

- 6) Do you require the 21 CFR Part 11 Module? Yes No
The xPONENT Software provides the user the opportunity to enable the 21CFR Part 11 module for regulatory compliance. This module maintains data integrity through the use of electronic signatures. The 21CFR Part 11 module contains the Basic Security module.

- 7) Do you require the Laboratory Information System (LIS) Module? Yes No
*This module enables **xPONENT 3.1** software to transmit and receive data from Laboratory Information Systems (LIS) and Laboratory Information Management Systems (LIMS) using an ASTM file format.*

- 8) Do you require the Automation Interface Module? Yes No
*The **xPONENT 3.1** software will facilitate the communication/integration between front-end automation devices and Luminex 100/200 systems. The front-end automation provider can build the required driver through an Application Programming Interface (API) which will be supplied with the Automation Interface Module.*

- 9) Do you require additional seat licenses (beyond the 3 that are standard)? Yes No
If so, how many additional seats do you require? _____
The maximum number of seat licenses per Luminex system is 15.

- 10) Do you require additional training (beyond the 1 web-based training course included)? Yes No
 Web-based training; number of seats _____ Instructor-led training; number of seats _____

Once you have completed the System Requirements Checklist and the Upgrade Checklist, provide it to your sales representative who will use it to create your quote. Please reference this quote number when you order your customized **xPONENT 3.1** upgrade kit.

We look forward to hearing from you!

Thank you,
Luminex Corporation