



# GETTING PROService

Please review the requirements below and complete the following questionnaire for PROService implementation.



Move healthcare forward.

## PROService network requirements and questionnaire

### Firewall

- > All inbound ports are closed
- > All outbound communications are restricted to only the Beckman Coulter domain
- > Required hospital firewalls with Beckman Coulter servers:
  - Outbound proservicedc.beckman.com (134.217.3.16) Port 443
  - Outbound proservicerds.beckman.com (134.217.3.17) Port 443
    - Allow OpenVPN on Port 443
    - Allow all subsequent established/connected traffic after the initial connection is established
- > Outbound www.beckmancoulter.com (134.217.3.37) Port 443
- > Instruments on the RAP Box subnet are blocked from exploring devices outside of the instrument subnet

### Instrument subnet

172.16.1.XXX or 192.168.1.XXX with a subnet mask of 255.255.255.0

### Optional Wi-Fi connection

- > Requires static network security authentication
- > Connects to any existing 2.4 or 5 GHz Wi-Fi access point, gateway or router
- > Supports Wi-Fi Protected Access\* (WPA2-PSK, WPA-PSK) and WEP data encryption

### Customer information

Facility name: \_\_\_\_\_

Ship-to address: \_\_\_\_\_

City: \_\_\_\_\_ State: \_\_\_\_\_ Zip code: \_\_\_\_\_ Country: \_\_\_\_\_

### Ship-to contact

Name: \_\_\_\_\_ Phone: \_\_\_\_\_

### IT contact

Name: \_\_\_\_\_

Phone: \_\_\_\_\_ Email: \_\_\_\_\_

### Laboratory contact

Name: \_\_\_\_\_

Phone: \_\_\_\_\_ Email: \_\_\_\_\_

### Network settings

#### Preferred network connection type:

Fast Ethernet hardwire

Is there a network line available to connect?

Yes  No  M jack #: \_\_\_\_\_

Optional wireless Wi-Fi  (Wireless connection requires static security authentication and does not support password rotation.)

Encryption type: WPA2-PSK  WPA-PSK

SSID: \_\_\_\_\_

Username: \_\_\_\_\_

Password: \_\_\_\_\_

**Network settings (cont.)**

**Does your firewall require a MAC address (select only one)?** Yes  No

Static IP  DHCP

IP address: \_\_\_\_\_ Subnet mask: \_\_\_\_\_

Default gateway: \_\_\_\_\_ DNS server 1: \_\_\_\_\_

DNS server 2 (optional): \_\_\_\_\_

**Is a proxy server used for Internet access (select only one)?** Yes  No

HTTP  SOCKS

IP address: \_\_\_\_\_

Port number: \_\_\_\_\_

Authenticating proxy (select only one)? Yes  No

Proxy username: \_\_\_\_\_ Proxy password: \_\_\_\_\_

**Is Port 443 available (select only one)?** Yes  No

**Connected instruments for PROService****Automation and middleware**

AutoMate 600  REMISOL Advance

**Chemistry**

DxC 600/800  DxC 700 AU  AU680  AU5800

**Hematology**

DxH 600/800  DxH SMS  LH500  LH750

**Immunoassay**

DxI 600/800  Access 2

**Life sciences**

AQUIOS  Biomek i-Series  Gallios  Navios

**Instance ID/Model/Serial number of instruments to be connected:**

To submit your completed questionnaire or request additional information, please email [proserviceinstalls@beckman.com](mailto:proserviceinstalls@beckman.com) or fax to (714) 223-4180.

REMISOL Advance is a trademark of Normand-Info SAS.

© 2017 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit [www.beckmancoulter.com/contact](http://www.beckmancoulter.com/contact)

FL-44712