



# Industrial Application Note

## 8011+ Web Server Application Setup Procedure

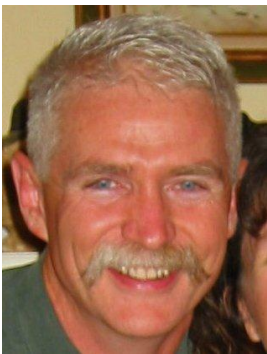
## **8011+ Web Server Application Setup Procedure**

1. Connect the 8011+ Ethernet port to your Network via a suitable Cat 5 Ethernet cable. If the cable is not connected first then the system will continue to scan the port until a cable is connected.
2. Power up the 8011+ and wait until initialization and self-tests are complete.
3. From the Home screen press **Setup**, then press **Instrument**, then press **Web Server**.
4. From the Web Server screen set the **Web Server**.
  - a. Option 1: Set to **Yes** by using the “1” button
  - b. Option 2: Web Server name – Create your personalized Web Server name (this will be the name of the 8011+ Web Server that will be indicated (seen) on the Network. The Default is: “**ABS-your 8011+ serial number**”. Example: “**ABS-1404554008**”)
  - c. Option 3: DHCP – Set this to **Yes** if you wish the IP Address to be automatically assigned by the network, or to **No** if you wish to manually set the **IP Address**, **Subnet Mask**, and **Gateway** values.
5. After this setup is complete, then press **Setup**, then **Home**, then press **Historical Data**, and then press **More**, and then press **Export all records**.
6. You will have 3 Options:
  1. Data to export - Defines data to be exported to the Web Server
    1. All data
    2. Filtered data only (this will have to be set up in the Historical data screen)
    3. Database file (Not Applicable for Web Server application)
  2. Output format
    1. TSV
    2. PDF
  3. Output Location
    1. Web Server (select Web Server)
    2. USB

7. When the options have been selected press **Export**.
8. The data file will now be created (see “**Preparing Export Data**” message at the bottom of the display screen). “**Export was successful**” message will appear on the screen when the activity is complete.
9. The Data file will now be available on the Web Server by the file name you created in Step 4.b.
10. To access the data go to a computer that is connected to the same Network and open a Web Browser like Internet Explorer.
11. **Type** in the URL path: “http://Your webserver name” (or the device IP address established in Step 4.c.) and then press **Enter**.
12. Your Web Server Data file should show up as a link and now you can click on it to access the data specified in step 6.a.

## **Author Biography**

Bill F. Bars is an Application Scientist for Beckman Coulter Life Sciences Company in Grants Pass, Oregon, USA. He has created and developed many of the Industrial Systems production processes and procedural tools for the Beckman Coulter / Hach Ultra Particle Counting Business Units products. These products include but are not limited to the: **HIAC PODS, 8011, 8011+, 8012, HRLD Sensors, PM4000, ROC, and the Calibration Fluids Lab.** He was a primary technical resource for the Hach Ultra Particle Counting ISO 17025 accreditation project which culminated in receiving their formal ISO 17025 Accreditation Certificate from A2LA. He received his Electronics Engineering degree from DeVry Institute of Technology in 1982. He has worked for the Beckman Coulter/Hach Ultra Companies for 17 years in a multitude of engineering capacities ranging from Metrology to Service Training and Industrial Application Support.



Beckman Coulter Life Sciences  
Particle Counting and Characterization  
481 California Ave  
Grants Pass, OR 97526 USA  
www.particle.com  
Information 1-800-866-7889, Ext 6195  
Email Bill Bars at: [bbars@beckman.com](mailto:bbars@beckman.com)