

# Automated, Accurate pH and Conductivity Measurements Using a Discrete Photometric Analyzer with an ECM Module

Mari Kiviluoma, Sari Hartikainen  
Thermo Fisher Scientific Oy, Analyzers & Automation, Vantaa, Finland

## Introduction

In this study pH and electrical conductivity (EC) were measured from water samples using an Electro Chemical Measurement (ECM) unit integrated into an automated discrete photometric analyzer. Integrated ECM unit is capable of simultaneously measure both pH and conductivity alongside the photometric testing. Accuracy of the results was verified by parallel testing with manual conductivity and pH meters.

Conductivity test has been developed to cover a measurement range from 20  $\mu\text{S}/\text{cm}$  to 112  $\text{mS}/\text{cm}$ . The pH test measures a pH range from 2 to 12. Results show very good precision, below 1.0 %, and good correlation to the manual methods with an  $R^2$  value ranging from 0.977 to 1.000.

## Methods

In the ECM unit, the conductivity measurement is performed via two electrodes. pH is measured using a two-electrode galvanic cell consisting of an indicator pH electrode and a reference electrode.

### Temperature compensation

Both conductivity and pH are measured at 37 °C. However, results can be reported in different temperatures, e.g., at 25 °C, because the discrete analyzer software has a robust system with which the sample result may be correlated to a reference analyzer result automatically.

Published correction factor values for conductivity measurements are available for a variety of sample types.

## Instruments

Thermo Scientific™ Gallery™ Plus analyzer with the ECM unit.

Thermo Scientific™ Orion™ 3 Star benchtop pH meter with Orion ROSS™ Combination Glass pH Electrode.

Radiometer™ CDM 80 Conductivity meter.

## Calibration

Calibration in the Gallery analyzer is performed using point-to-point calibration type. With a point-to-point calibration separate linear equations are generated between each set of two calibrators. For example if the calibration is done using five calibration levels, four separate equations are generated for the four linear segments.

FIGURE 1. Conductivity calibration.

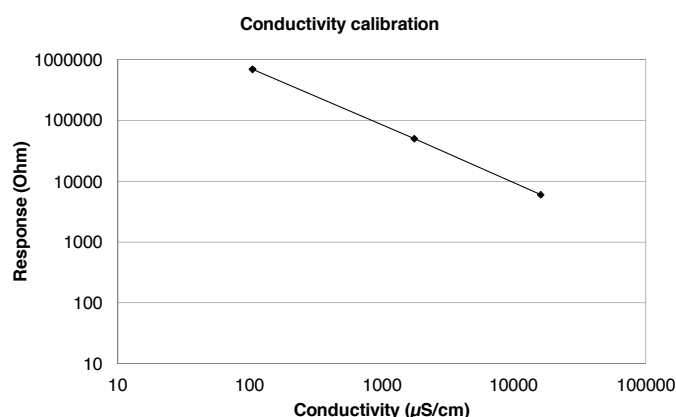


FIGURE 2. pH calibration.

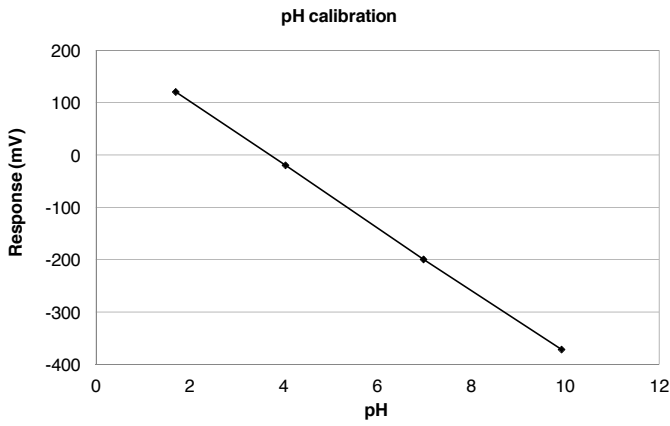
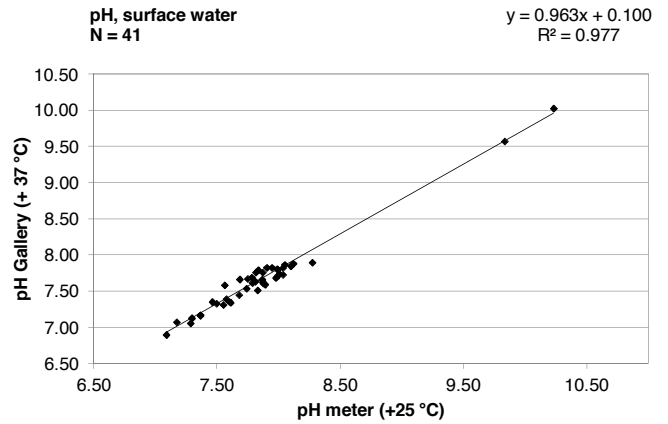


FIGURE 4. Method comparison of surface water samples.



## Results

All measurements were done simultaneously with the Gallery Plus discrete analyzer and the manual pH or conductivity meter. Figures from 3 to 6 shows the method comparison results. Temperature compensation is not included in the graphs, but correction factors and biases for different matrices are calculated to the Table 1.

FIGURE 3. Method comparison of natural water samples containing surface and ground water samples. Samples were measured at +37°C using Gallery temperature compensation with the correction factor of 0.793 (from literature).

FIGURE 5. Method comparison of ground and well water samples.

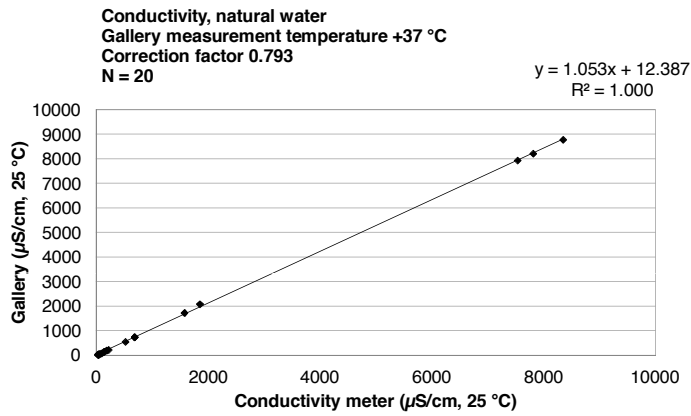
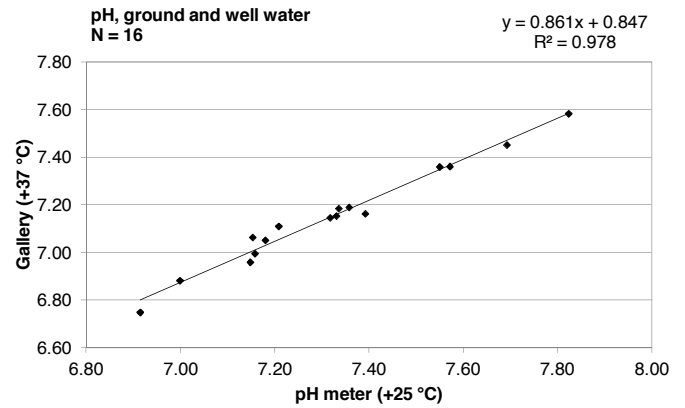
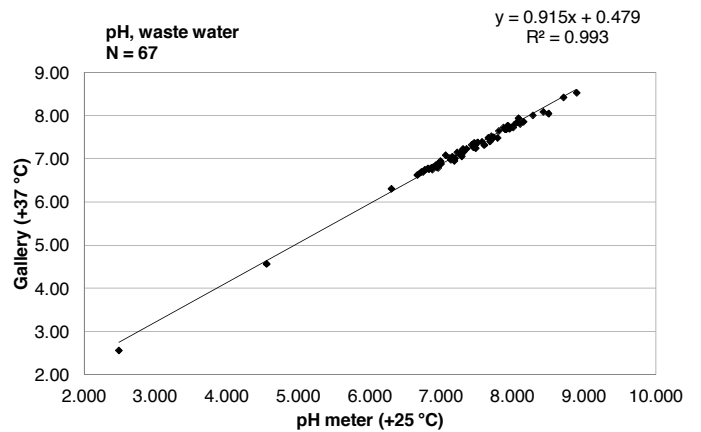


FIGURE 6. Method comparison of waste water samples containing influents, effluents and industrial waste waters.



**TABLE 1.** Calculated correction factors and biases for different matrices based on method comparison results.  $C_{factor}$  is the inverse of the slope and  $C_{bias}$  is the unmodified y-intercept.

	$C_{factor}$	$C_{bias}$
Surface water	1.038	0.100
Ground and well water	1.161	0.847
Waste water	1.093	0.479

**TABLE 2.** Precision results of conductivity and pH.

	Natural water			
	Conductivity		pH	
	N	20	N	20
	Mean ( $\mu\text{S/cm}$ )	194.13	Mean	7.80
	SD	CV %	SD	CV %
<b>Within Run</b>	0.709	0.37 %	0.061	0.78 %
<b>Between Run</b>	1.410	0.73 %	0.007	0.09 %
<b>Total</b>	1.578	0.81 %	0.062	0.79 %

## Conclusion

The automated pH and conductivity methods correlate well with the manual methods. Although measurement temperature in the ECM unit is higher, the temperature compensation can be done automatically with the software and the results can be reported at +25°C.

## References

1. Laxen, D.P.H., A specific conductance method for quality control in water analyses. In: Water Research, Vol 11, 1977, pp 91-94.
2. SFS-EN 27888:1985, Suomen Standardisoimisliitto SFS. Water quality. Determination of electrical conductivity.

[www.thermoscientific.com](http://www.thermoscientific.com)

©2016 Thermo Fisher Scientific Inc. All rights reserved. Radiometer is the trademark of Danaher Corporation. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.



Thermo Fisher Scientific,  
Sunnyvale, CA USA is  
ISO 9001 Certified.

**Africa** +43 1 333 50 34 0  
**Australia** +61 3 9757 4300  
**Austria** +43 810 282 206  
**Belgium** +32 53 73 42 41  
**Brazil** +55 11 2730 3006  
**Canada** +1 800 530 8447  
**China** 800 810 5118 (free call domestic)  
 400 650 5118

**Denmark** +45 70 23 62 60  
**Europe-Other** +43 1 333 50 34 0  
**Finland** +358 10 3292 200  
**France** +33 1 60 92 48 00  
**Germany** +49 6103 408 1014  
**India** +91 22 6742 9494  
**Italy** +39 02 950 591

**Japan** +81 6 6885 1213  
**Korea** +82 2 3420 8600  
**Latin America** +1 561 688 8700  
**Middle East** +43 1 333 50 34 0  
**Netherlands** +31 76 579 55 55  
**New Zealand** +64 9 980 6700  
**Norway** +46 8 556 468 00

**Russia/CIS** +43 1 333 50 34 0  
**Singapore** +65 6289 1190  
**Sweden** +46 8 556 468 00  
**Switzerland** +41 61 716 77 00  
**Taiwan** +886 2 8751 6655  
**UK/Ireland** +44 1442 233555  
**USA** +1 800 532 4752

PN71982-EN 0216S

**Thermo**  
SCIENTIFIC

A Thermo Fisher Scientific Brand