

Measuring pH in Low Ionic Strength Solutions

Water Analysis Instruments, Thermo Fisher Scientific

Key Words

pH, low ionic strength, pure water, electrodes, solutions, ROSS, ROSS Ultra, water conductivity, pHISA adjuster, pure water pH test kit.

Goal

The following application note describes the challenges and recommendations when measuring pH in low ionic strength solutions.

Introduction

pH measurements are typically made in solutions which contain relatively large amounts of acid or base, or which contain substantial amounts of dissolved salts. Under these conditions, conventional pH electrodes make measurements with speed and precision.

There is a strong interest in making pH measurements in “pure water” (for example, waters in which the total amount of acid or base is very small, and in which there is a low level of dissolved salts). The terms “pure water” and “low ionic strength” can be used interchangeably. Samples which may fall into this category include:

- Distilled waters
- Deionized waters
- Some process waters
- Well waters
- Some surface waters
- Treated effluent water
- Boiler feed water

Measuring pH in these pure water samples can be difficult. Although electrodes respond quickly in buffers, their response in pure water can be unsatisfactory – slow, noisy, drifting, non-reproducible, and inaccurate.

The Problem

These commonly observed problems can be attributed to the low conductivity of the sample, differences between low ionic strength solutions and normal ionic strength buffers, changes in the liquid junction potential, and absorption of carbon dioxide.



Because pure water samples are poor conductors, the solution will often behave like “antenna,” and the electrode response can be noisy.

Standardization of an electrode in a high ionic strength buffer will increase the time required for stabilization when going to a low ionic strength solution. In addition, the possibility of sample contamination will be increased. To ensure the most precise measurements, buffers and samples should close in ionic strength.

When any two solutions come in contact, diffusion occurs until equilibrium is reached. Since ions have different mobilities and diffuse at different rates, a charge imbalance occurs causing a junction potential. A junction potential occurs when the reference electrode fill solution meets the sample.

This will be large if there is a large difference in composition between the fill solution and the sample. It is important that the junction potential be constant during measurement. If the two solutions are quite different, normal fluctuations in the boundary layer will produce noise. Constant, reproducible junction potentials are achieved by measuring in samples and standards with ionic strengths similar to the fill solution.

Since pure waters contain little dissolved material, their buffering capacity is small. Absorption of carbon dioxide from the atmosphere will cause a slow change in the pH, which is observed as a drift in pH reading and a different pH from the original sample. Samples which are not previously saturated with carbon dioxide must be handled with care.

Conventional Approaches

The most widely accepted approaches to solving these problems involve the use of low resistance glass pH electrodes or the use of a reference electrode with a fast, continuous leak rate. When placed in a pure water sample, these electrodes show improved time response and stability, due to dissolution of the low resistance glass into the low ionic strength solution and the non-quantitative addition of a salt solution from the reference into the sample respectively.

Both techniques raise the conductivity, but, at the same time, may change the sample pH at the electrode surface. Response is improved, but an error is added to the measurement, which depends upon the amount of dissolved material. Increased sample conductivity is desirable; however, non-reproducible alteration of the pH reading is not.

To purchase an Orion ROSS electrode or the Orion Pure Water pH Test Kit, please contact your local equipment distributor and reference the part numbers listed below:

Product	Description	Part Number
Electrode	Thermo Scientific Orion ROSS Ultra pH Electrode	8102BNUWP
	Thermo Scientific™ Orion™ ROSS Ultra™ Glass Triode™ pH/ATC Combination Electrode	8302BNUMD
	Thermo Scientific Orion ROSS Ultra Low Maintenance pH/ATC Triode Combination Electrode	8107BNUMD/8107UWMMD
	Thermo Scientific Orion ROSS Ultra Refillable pH/ATC Triode Combination Electrode	8157BNUMD/8157UWMMD
Accessories	Thermo Scientific Orion Pure Water pH Test Kit	700001

Recommended Electrodes and Solution

When measuring pH in low ionic strength solutions, the following equipment is recommended to overcome the challenges presented by the conventional approaches:

- Thermo Scientific™ Orion™ ROSS™ electrode or Thermo Scientific™ Orion™ ROSS Ultra™ electrode
- Thermo Scientific™ Orion™ Pure Water™ pH Test Kit

The kit includes the Thermo Scientific™ Orion™ pHISA™, a pure water pH additive and a special set of buffers containing the same background of pHISA adjustor. Adding pHISA adjustor to samples increases the ionic strength, thus reducing noise and improving response time. The shift in pH caused by the addition of pHISA adjustor is minimal, between 0.005 and 0.01 pH units. Since the same amount of pHISA adjustor is added to the buffers and samples, the net effect on the pH is negligible.

Conclusion

Calibration is performed using special dilute buffers with pHISA adjustor already added. Measuring with samples and buffers of the same ionic strength increases accuracy, precision, and response time. Contamination due to carryover from higher ionic strength buffers is also minimized.

Errors in pH measurement due to liquid junction potential variations are minimized by using buffers and samples at similar ionic strength. Addition of pHISA adjustor to both the buffers and samples achieves this condition.

Junction potentials will vary depending on the style of junction and choice of electrode fill solution. A good quality pH electrode will provide better response when compared against the universal standard, the hydrogen electrode. To optimize the measurement, we recommend using the same Orion ROSS electrode.

thermoscientific.com/water

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries.

Water Analysis Instruments

North America
Toll Free: 1-800-225-1480
Tel: 1-978-232-6000
info.water@thermo.com

Netherlands
Tel: (31) 020-4936270
info.water.uk@thermo.com

China
Tel: (86) 21-68654588
wai.asia@thermofisher.com

India
Tel: (91) 22-4157-8800
wai.asia@thermofisher.com

Singapore
Tel: (65) 6778-6876
wai.asia@thermofisher.com

Japan
Tel: (81) 045-453-9175
wai.asia@thermofisher.com

Australia
Tel: (613) 9757-4300
in Australia (1300) 735-295
InfoWaterAU@thermofisher.com

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand