

Orion AquaPro Series Combination pH Electrode Instruction Sheet

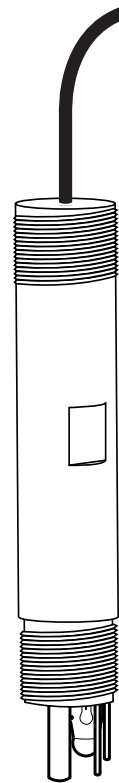
AQUA10 pH Electrode

The AQUA10 is an industrial combination pH electrode designed for accurate pH measurements in the harshest environments and samples. The AQUA10 is a double junction Ag/AgCl non-refillable pH electrode with automatic temperature compensation and solution ground. Three stainless steel bulb guards protect the pH sensing bulb from particulates or solids in a flowing sample. The electrode has $\frac{3}{4}$ " NPT threads for mounting within pipe systems or at the end of a piece of conduit for submersion applications. The AQUA10 pH electrode is compatible for use with the Orion 2001 pH Monitor.

For Best Results:

- Inspect the electrode for scratches, cracks, salts or other deposits on the sensing element and reference junction. Rinse off any precipitates with distilled water.
- To attach the electrode to the monitor, wire the stripped, unterminated ends to the monitor. Refer to the monitor instruction manual.
- The first time a new electrode is installed, condition the electrode in pH 7 buffer (Orion 200107) for 1 hour.
- Put the electrode on-line and allow the electrode to equilibrate in the sample for a few hours prior to calibration.
- Always use fresh buffers when calibrating. Use Orion precision buffers: pH 4 (Orion 200104), pH 7 (Orion 200107) or pH 10 (Orion 200110). Choose buffers that bracket the expected sample pH. Check electrode slope by performing a two-point buffer calibration. Slope should be within 92 to 102 %. For detailed calibration and temperature compensation procedures, consult the Orion 2001 pH Monitor instruction manual.
- To maximize the life of your pH electrode, use Orion pH Electrode Cleaning Kits. After cleaning, soak the electrode in pH 7 buffer (200107) for at least 1 hour. Recalibrate before returning to on-line measurement.

Note: see reverse side for specifications and ordering information



Specifications:

pH Range:	0 to 14
Temperature Range:	-5 to 60 °C
Pressure Range:	0 to 100 psi
Drift:	≤ 0.5 pH units for 30 days
Reproducibility:	± 0.01 pH units
Slope:	92 to 102 % (54.4 to 60.3 mV/dec)
ATC Resistance:	30 kΩ
Reference System:	<i>CONTACT</i> Reference System
Reference Electrolyte:	non-refillable, <i>CONTACT</i> polymer gel
AQUA10 Diameter:	30 mm
AQUA10 Length:	170 mm
Cable:	coaxial, 5 meters

Ordering Information

Orion	Description
AQUA10	Industrial Combination pH Electrode.
900021	pH Cleaning Solution A, for removing protein contaminants. Four 1 oz. bottles
900022	pH Cleaning Solution B, for general cleaning and bacterial contaminants. Four 2 oz. bottles
900023	pH Cleaning Solution C, for general cleaning. Four 1 oz. bottles
900024	pH Cleaning Solution D, for oil and grease contaminants. Four 2 oz. bottles

Environmental Instruments Water Analysis

For updated contact information,
visit www.thermo.com

North America
166 Cummings Center
Beverly, MA 01915 USA
Tel: 978-232-6000
Dom. Fax: 978-232-6015
Int'l. Fax: 978-232-6031

Europe
12-16 Sedgeway Business Park
Witchford, Cambridgeshire
England, CB6 2HY
Tel: 44-1353-666111
Fax: 44-1353-666001

Far East
Room 904, Federal Building
369 Lockhart Road
Wanchai, Hong Kong
Tel: 852-2836-0981
Fax: 852-2834-5160

Customer Support
Toll Free: 800-225-1480
www.thermo.com
Dom. e-mail: domcs1@thermoorion.com
Int'l. e-mail: intcs1@thermoorion.com