

Programming Data for Spectroquant® Test Kits

Calibration Information: wavelength, slope and blank value

For swift, secure analysis, there's no better choice than Spectroquant® test kits. Consisting of validated, standard-compliant reagents, the kits are pre-programmed for use with Spectroquant® instruments to ensure rapid, reliable results. As a result of their excellent quality, most of our kits conform to international standards, allowing you to test with absolute confidence.

Spectroquant® Reagent Tests contain all reagents for your photometric analyses and can be used with any existing cells of various pathlengths (10 - 100-mm rectangular cells or 16-mm / 1-inch round cells). The measurement range and programming data depend on the size of the cells (which must be acquired separately). In the table below, these Reagent Tests are abbreviated as **RT**.

Spectroquant® Cell Tests are ready-to-use 16-mm (OD) round cells pre-filled with reagents; no additional cell is needed. They are even more easy-to-use and convenient. In the table below, these Cell Tests are abbreviated as **CT**.

All test kits can be used seamlessly with **Spectroquant® Photometers**. Calibration data are pre-programmed, and the workflow is designed to enable fast and reliable results with additional features for convenience. For example, the **Spectroquant® Prove** photometer series can read Live ID barcodes that provide batch-specific information such as the shelf-life. It also automatically recognizes the inserted cell and adapts the measurement range to display the correct result without any action needed from the user. The system also includes ready-to-use reference materials and is designed to comply with highest requirements in terms of analytical quality assurance. For further details, please visit www.sigmaaldrich.com/photometry.



If you already have a photometer in your lab and want to experience the outstanding quality and reliability of the Spectroquant® test kit series, you may use the programming data provided in this document. Please make sure that the instruments can (a) measure at the listed wavelength and (b) hold the respective cell sizes as there can be different models available (especially relevant for the round cell dimensions) and if e.g., an adapter is needed.

The calibration data were obtained on different reference photometers from different manufacturers at a certain point in time. Please note that we cannot guarantee the quality of measurement results in this case, as we neither monitor potential changes in the instruments from other manufacturers nor control the performance of your photometer in terms of wavelength precision, repeatability, and linearity of absorbance. In any case we recommend to validate each new method with standards after programming it on your device. But if your instrument is high-quality, and is regularly subjected to the appropriate quality assurance (e.g. with Supelco® reference standards), Spectroquant® test kits can provide a unique experience in terms of precision, convenience, and compliance.

When programming a method you will also need to measure against a blank in most cases. The blanks the below methods are calibrated for can be:

- Dist. Water (DW) means distilled water without reagents and poured into the same type of cell used for the measurement.
- Own blank (OB) means distilled water and all reagents in the same quantity as used for the measuring sample preparation (can also be called reagent blank). Separately prepared in standard test tubes and poured into the same type of cell used for the measurement.
- Sample blank (SB) means to measure the sample material as a blank (turbid samples have to be filtered), without reagents and poured into the same type of cell used for the measurement.

If the calibration function of the respective test kit is not linear, you will find the remark A) in the factor column, and the respective calibration information is given from page 18.

If you need more information, please contact our technical service at www.sigmaaldrich.com, or your trusted dealer.

Programming Data

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|-------------------------|-----------------------------|-----------|---------------------|-------------|--------|------------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Acid Capacity to pH 4.3 | 1.01758 | CT | 0.40 | 8.00 | mmol/l | OH ⁻ | 16 mm | A) | 605 | DW |
| | | | 0.40 | 8.00 | mmol/l | OH ⁻ | 10 mm | | | |
| | | | 20 | 400 | mg/l | CaCO ₃ | 16 mm | | | |
| | | | 20 | 400 | mg/l | CaCO ₃ | 10 mm | | | |
| Acid Capacity to pH 4.3 | 1.01762 | CT | replaced by 1.01758 | | | | | | | |
| Alcohol | Test deleted | | | | | | | | | |
| Alkalinity | see Acid capacity to pH 4.3 | | | | | | | | | |
| Aluminium | 1.14825 | RT | 0.10 | 1.20 | mg/l | Al | 10 mm | A) | 550 | DW |
| | | | 0.020 | 0.200 | mg/l | Al | 50 mm | | | |
| | | | 0.039 | 0.472 | mg/l | Al | 1 inch | | | |
| Aluminium | 1.00594 | CT | 0.02 | 0.50 | mg/l | Al | 16 mm | A) | 545 | DW |
| | | | 0.02 | 0.50 | mg/l | Al | 10 mm | | | |
| Ammonium | 1.14739 | CT | 0.010 | 2.000 | mg/l | NH ₄ -N | 16 mm | 0.877 | 690 | OB |
| | | | 0.010 | 2.000 | mg/l | NH ₄ -N | 10 mm | 1.18 | | |
| | | | 0.01 | 2.58 | mg/l | NH ₄ ⁺ | 16 mm | 1.13 | | |
| | | | 0.01 | 2.58 | mg/l | NH ₄ ⁺ | 10 mm | 1.53 | | |
| | | | 0.010 | 2.000 | mg/l | NH ₃ -N | 16 mm | 0.877 | | |
| | | | 0.010 | 2.000 | mg/l | NH ₃ -N | 10 mm | 1.18 | | |
| | | | 0.01 | 2.43 | mg/l | NH ₃ | 16 mm | 1.065 | | |
| 0.01 | 2.43 | mg/l | NH ₃ | 10 mm | 1.433 | | | | | |
| Ammonium | 1.14752 | RT | 0.05 | 3.00 | mg/l | NH ₄ -N | 10 mm | 1.23 | 690 | OB |
| | | | 0.03 | 1.50 | mg/l | NH ₄ -N | 20 mm | 0.615 | | |
| | | | 0.010 | 0.500 | mg/l | NH ₄ -N | 50 mm | 0.246 | | |
| | | | 0.02 | 1.18 | mg/l | NH ₄ -N | 1 inch | 0.483 | | |
| | | | 0.06 | 3.86 | mg/l | NH ₄ ⁺ | 10 mm | 1.58 | | |
| | | | 0.04 | 1.93 | mg/l | NH ₄ ⁺ | 20 mm | 0.792 | | |
| | | | 0.013 | 0.644 | mg/l | NH ₄ ⁺ | 50 mm | 0.316 | | |
| | | | 0.024 | 1.520 | mg/l | NH ₄ ⁺ | 1 inch | 0.622 | | |
| | | | 0.05 | 3.00 | mg/l | NH ₃ -N | 10 mm | 1.23 | | |
| | | | 0.03 | 1.50 | mg/l | NH ₃ -N | 20 mm | 0.615 | | |
| | | | 0.010 | 0.500 | mg/l | NH ₃ -N | 50 mm | 0.246 | | |
| | | | 0.02 | 1.18 | mg/l | NH ₃ -N | 1 inch | 0.483 | | |
| | | | 0.06 | 3.65 | mg/l | NH ₃ | 10 mm | 1.494 | | |
| | | | 0.04 | 1.82 | mg/l | NH ₃ | 20 mm | 0.747 | | |
| | | | 0.016 | 0.608 | mg/l | NH ₃ | 50 mm | 0.384 | | |
| 0.03 | 1.44 | mg/l | NH ₃ | 1 inch | 0.588 | | | | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

A) For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|-----------------|------------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Ammonium | 1.14558 | CT | 0.20 | 8.00 | mg/l | NH ₄ -N | 16 mm | 4.26 | 690 | OB |
| | | | 0.20 | 8.00 | mg/l | NH ₄ -N | 10 mm | 5.75 | | |
| | | | 0.26 | 10.30 | mg/l | NH ₄ ⁺ | 16 mm | 5.49 | | |
| | | | 0.26 | 10.30 | mg/l | NH ₄ ⁺ | 10 mm | 7.41 | | |
| | | | 0.20 | 8.00 | mg/l | NH ₃ -N | 16 mm | 4.26 | | |
| | | | 0.20 | 8.00 | mg/l | NH ₃ -N | 10 mm | 5.75 | | |
| | | | 0.24 | 9.73 | mg/l | NH ₃ | 16 mm | 5.17 | | |
| | | | 0.24 | 9.73 | mg/l | NH ₃ | 10 mm | 6.98 | | |
| Ammonium | 1.14544 | CT | 0.5 | 16.0 | mg/l | NH ₄ -N | 16 mm | 7.84 | 690 | OB |
| | | | 0.5 | 16.0 | mg/l | NH ₄ -N | 10 mm | 10.6 | | |
| | | | 0.6 | 20.6 | mg/l | NH ₄ ⁺ | 16 mm | 10.1 | | |
| | | | 0.6 | 20.6 | mg/l | NH ₄ ⁺ | 10 mm | 13.6 | | |
| | | | 0.5 | 16.0 | mg/l | NH ₃ -N | 16 mm | 7.84 | | |
| | | | 0.5 | 16.0 | mg/l | NH ₃ -N | 10 mm | 10.6 | | |
| | | | 0.6 | 19.5 | mg/l | NH ₃ | 16 mm | 9.52 | | |
| | | | 0.6 | 19.5 | mg/l | NH ₃ | 10 mm | 12.87 | | |
| Ammonium | 1.00683 | RT | 2.0 | 75.0 | mg/l | NH ₄ -N | 10 mm | 27.8 | 690 | OB |
| | | | 5 | 150 | mg/l | NH ₄ -N | 10 mm | 55.6 | | |
| | | | 2.6 | 96.6 | mg/l | NH ₄ -N | 10 mm | 35.8 | | |
| | | | 6 | 193 | mg/l | NH ₄ ⁺ | 10 mm | 71.5 | | |
| | | | 2.0 | 75.0 | mg/l | NH ₃ -N | 10 mm | 27.8 | | |
| | | | 5 | 150 | mg/l | NH ₃ -N | 10 mm | 55.6 | | |
| | | | 2.4 | 91.2 | mg/l | NH ₃ | 10 mm | 33.8 | | |
| | | | 6 | 182 | mg/l | NH ₃ | 10 mm | 67.5 | | |
| Ammonium | 1.14559 | CT | 4.0 | 80.0 | mg/l | NH ₄ -N | 16 mm | 36.4 | 690 | OB |
| | | | 4.0 | 80.0 | mg/l | NH ₄ -N | 10 mm | 49.1 | | |
| | | | 5.2 | 103.0 | mg/l | NH ₄ ⁺ | 16 mm | 46.8 | | |
| | | | 5,2 | 103.0 | mg/l | NH ₄ ⁺ | 10 mm | 63.3 | | |
| | | | 4.0 | 80.0 | mg/l | NH ₃ -N | 16 mm | 36.4 | | |
| | | | 4.0 | 80.0 | mg/l | NH ₃ -N | 10 mm | 49.1 | | |
| | | | 4.9 | 97.3 | mg/l | NH ₃ | 16 mm | 44.2 | | |
| | | | 97.3 | mg/l | NH ₃ | 10 mm | 59.6 | | | |
| AOX | 1.00675 | RT | 0.05 | 2.50 | mg/l | AOX | 16 mm | 4.00 | 445 | OB |
| | | | 0.05 | 2.50 | mg/l | AOX | 10 mm | 5.40 | | |
| Antimony | C) | | 0.10 | 8.00 | mg/l | Sb | 10 mm | 3.45 | 620 | OB |
| Arsenic | 1.01747 | RT | 0.005 | 0.100 | mg/l | As | 10 mm | 0.1138 | 525 | DW |
| | | | 0.001 | 0.020 | mg/l | As | 20 mm | 0.0568 | | |
| | | | 0.002 | 0.039 | mg/l | As | 1 inch | 0.0447 | | |
| BOD | 1.00687 | CT | 0.5 | 3000 | mg/l | BOD | 16 mm | 13.5 | 500 | DW |
| | | | 0.5 | 3000 | mg/l | BOD | 10 mm | 18.2 | | |
| Boron | 1.14839 | RT | 0.050 | 0.800 | mg/l | B | 10 mm | 0.328 | 565 | OB |
| Boron | 1.00826 | CT | 0.05 | 2.00 | mg/l | B | 16 mm | 1.38 | 405 | OB |
| | | | 0.05 | 2.00 | mg/l | B | 10 mm | 1.86 | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaldrich.com.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|---------------------------------|-----------------------|-----------|-----------------|-------------|------------|-------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Bromine | E) | RT | 0.10 | 10.00 | mg/l | Br₂ | 10 mm | 7.52 | 550 | DW |
| | | | 0.05 | 5.00 | mg/l | Br₂ | 20 mm | 3.76 | | |
| | | | 0.020 | 2.000 | mg/l | Br₂ | 50 mm | 1.50 | | |
| | | | 0.039 | 3.937 | mg/l | Br₂ | 1 inch | 2.960 | | |
| Cadmium | 1.01745 | RT | 0.010 | 0.500 | mg/l | Cd | 10 mm | 0.680 | 525 | OB |
| | | | 0.005 | 0.250 | mg/l | Cd | 20 mm | 0.340 | | |
| | | | 0.002 | 0.100 | mg/l | Cd | 50 mm | 0.136 | | |
| | | | 0.005 | 0.197 | mg/l | Cd | 1 inch | 0.268 | | |
| Cadmium | 1.14834 | CT | 0.025 | 1.000 | mg/l | Cd | 16 mm | 0.595 | 550 | OB |
| | | | 0.025 | 1.000 | mg/l | Cd | 10 mm | 0.804 | | |
| Calcium | 1.00049 | RT | 0.20 | 4.00 | mg/l | Ca | 10 mm | D) | 565 | OB |
| Calcium | 1.14815 | RT | 10 | 160 | mg/l | Ca | 10 mm | 137 | 550 | OB |
| | | | 25 | 400 | mg/l | CaCO₃ | 10 mm | 342 | | |
| | | | 14 | 224 | mg/l | CaO | 10 mm | 192 | | |
| | | | 5 | 80 | mg/l | Ca | 20 mm | 68.5 | | |
| | | | 12 | 200 | mg/l | CaCO₃ | 20 mm | 171 | | |
| | | | 7 | 112 | mg/l | CaO | 20 mm | 95.9 | | |
| | | | 1.0 | 15.0 | mg/l | Ca | 10 mm | 29.4 | | |
| | | | 2.5 | 37.5 | mg/l | CaCO₃ | 10 mm | 73.5 | | |
| Calcium | 1.00858 | CT | 10 | 250 | mg/l | Ca | 16 mm | A) | 565 | DW |
| | | | 25 | 625 | mg/l | CaCO₃ | 16 mm | | | |
| | | | 14 | 350 | mg/l | CaO | 16 mm | | | |
| | | | 10 | 250 | mg/l | Ca | 10 mm | | | |
| | | | 25 | 625 | mg/l | CaCO₃ | 10 mm | | | |
| | | | 350 | mg/l | CaO | 10 mm | | | | |
| Carbohydrazide | see Oxygen Scavengers | | | | | | | | | |
| Chloride | 1.01807 | RT | 0.10 | 5.00 | mg/l | Cl⁻ | 50 mm | A) | 500 | OB |
| Chloride | 1.01804 | CT | 0.5 | 15.0 | mg/l | Cl⁻ | 16 mm | A) | 445 | DW |
| | | | 0.5 | 15.0 | mg/l | Cl⁻ | 10 mm | | | |
| Chloride | 1.14897 | RT | 10 | 250 | mg/l | Cl⁻ | 10 mm | 107 | 500 | OB |
| | | | 2.5 | 125 | mg/l | Cl⁻ | 10 mm | 28.2 | | |
| Chloride | 1.14730 | CT | 5 | 125 | mg/l | Cl⁻ | 16 mm | 114 | 525 | OB |
| | | | 5 | 125 | mg/l | Cl⁻ | 10 mm | 153 | | |
| Chlorine (free chlorine) | 1.00598 | RT | 0.05 | 6.00 | mg/l | Cl₂ | 10 mm | 3.33 | 550 | DW |
| | | | 0.02 | 3.00 | mg/l | Cl₂ | 20 mm | 1.67 | | |
| | | | 0.010 | 1.000 | mg/l | Cl₂ | 50 mm | 0.667 | | |
| | | | 0.02 | 3.00 | mg/l | Cl₂ | 1 inch | 1.312 | | |
| Chlorine (free chlorine) | 1.00595 | CT | 0.03 | 6.00 | mg/l | Cl₂ | 16 mm | 2.47 | 550 | DW |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

A) For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.

D) Instrument-specific calibration needed. See Appendix page 19.

E) Can be determined with Chlorine test, 1.00598 (see corresponding application notes on www.sigmaaldrich.com).

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|---|-----------|-----------|---|-------------|------|--------------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Chlorine (total chlorine) | 1.00602 | RT | 0.05 | 6.00 | mg/l | Cl ₂ | 10 mm | 3.33 | 550 | DW |
| | | | 0.02 | 3.00 | mg/l | Cl ₂ | 20 mm | 1.67 | | |
| | | | 0.010 | 1.000 | mg/l | Cl ₂ | 50 mm | 0.667 | | |
| | | | 0.02 | 2.36 | mg/l | Cl ₂ | 1 inch | 1.312 | | |
| Chlorine (free + total chlorine) | 1.00597 | CT | 0.03 | 6.00 | mg/l | Cl ₂ | 16 mm | 2.47 | 550 | DW |
| Chlorine (free + total chlorine) | 1.00599 | RT | 0.05 | 6.00 | mg/l | Cl ₂ | 10 mm | 3.33 | 550 | DW |
| | | | 0.02 | 3.00 | mg/l | Cl ₂ | 20 mm | 1.67 | | |
| | | | 0.010 | 1.000 | mg/l | Cl ₂ | 50 mm | 0.667 | | |
| | | | 0.02 | 2.36 | mg/l | Cl ₂ | 1 inch | 1.312 | | |
| Chlorine (liquid) (free + total chlorine) | 1.00086 | RT | 0.03 | 6.00 | mg/l | Cl ₂ | 16 mm | 2.47 | 550 | DW |
| | 1.00087 | | 0.010 | 1.000 | mg/l | Cl ₂ | 50 mm | 0.667 | | |
| | 1.00088 | | 0.010 | 1.000 | mg/l | Cl ₂ | 50 mm | 0.667 | | |
| | 1.00089 | | 0.010 | 1.000 | mg/l | Cl ₂ | 50 mm | 0.667 | | |
| Chlorine | 1.14828 | RT | replaced by 1.00599 or 1.00598 or 1.00602 | | | | | | | |
| Chlorine | 1.19254 | RT | 0 | 2.00 | mg/l | Cl ₂ | 10 mm | B) | 520 or 530 | OB |
| | | | 0 | 2.00 | mg/l | Cl ₂ | 1 inch | B) | 520 or 530 | |
| Chlorine Dioxide | 1.00608 | RT | 0.10 | 10.00 | mg/l | ClO ₂ | 10 mm | 6.33 | 550 | DW |
| | | | 0.05 | 5.00 | mg/l | ClO ₂ | 20 mm | 3.16 | | |
| | | | 0.020 | 2.000 | mg/l | ClO ₂ | 50 mm | 1.27 | | |
| | | | 0.04 | 3.94 | mg/l | ClO ₂ | 1 inch | 2.49 | | |
| Chlorine Dioxide | 1.14732 | | replaced by 1.00608 | | | | | | | |
| Chromate | 1.14758 | RT | 0.05 | 3.00 | mg/l | Cr | 10 mm | 1.30 | 550 | DW |
| | | | 0.03 | 1.50 | mg/l | Cr | 20 mm | 0.650 | | |
| | | | 0.010 | 0.600 | mg/l | Cr | 50 mm | 0.260 | | |
| | | | 0.020 | 1.181 | mg/l | Cr | 1 inch | 0.511 | | |
| | | | 0.11 | 6.69 | mg/l | CrO ₄ ²⁻ | 10 mm | 2.90 | | |
| | | | 0.07 | 3.35 | mg/l | CrO ₄ ²⁻ | 20 mm | 1.45 | | |
| | | | 0.02 | 1.34 | mg/l | CrO ₄ ²⁻ | 50 mm | 0.579 | | |
| Chromate | 1.14552 | CT | 0.05 | 2.00 | mg/l | Cr | 16 mm | 0.971 | 550 | DW |
| | | | 0.05 | 2.00 | mg/l | Cr | 10 mm | 1.31 | | |
| | | | 0.11 | 4.46 | mg/l | CrO ₄ ²⁻ | 16 mm | 2.17 | | |
| | | | 0.11 | 4.46 | mg/l | CrO ₄ ²⁻ | 10 mm | 2.92 | | |
| Chromium Bath | C) | | 20 | 400 | g/l | CrO ₃ | 10 mm | 556 | 445 | DW |
| | | | 10 | 200 | g/l | CrO ₃ | 20 mm | 278 | | |
| | | | 4.0 | 80.0 | g/l | CrO ₃ | 50 mm | 111 | | |
| Cobalt | 1.17244 | CT | 0.05 | 2.00 | mg/l | Co | 16 mm | 3.715 | 500 | OB |
| | | | 0.05 | 2.00 | mg/l | Co | 10 mm | 5.02 | | |
| COD | 1.14560 | CT | 4.0 | 40.0 | mg/l | COD | 16 mm | -41.7 | 340 | OB |
| | | | 4.0 | 40.0 | mg/l | COD | 10 mm | -56.3 | | |
| COD | 1.01796 | CT | 5.0 | 80.0 | mg/l | COD | 16 mm | -40.7 | 340 | OB |
| | | | 5.0 | 80.0 | mg/l | COD | 10 mm | -54.9 | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

B) This test can (also) be used with Hach® factory-programmed instrument calibrations. No calibration factor/table is necessary. For the corresponding Hach program number see Appendix page 18.

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaaldrich.com.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|------------------------|------------------------|-----------|-----------------|-------------|-----------------|---------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| COD | 1.14540 | CT | 10 | 150 | mg/l | COD | 16 mm | -210 | 445 | OB |
| | | | 10 | 150 | mg/l | COD | 10 mm | -284 | | |
| COD | 1.14895 | CT | 15 | 300 | mg/l | COD | 16 mm | -222 | 445 | OB |
| | | | 15 | 300 | mg/l | COD | 10 mm | -300 | | |
| COD | 1.14690 | CT | 50 | 500 | mg/l | COD | 16 mm | -397 | 445 | OB |
| | | | 50 | 500 | mg/l | COD | 10 mm | -536 | | |
| COD | 1.14541 | CT | 25 | 1500 | mg/l | COD | 16 mm | 1667 | 605 | OB |
| | | | 25 | 1500 | mg/l | COD | 10 mm | 2249 | | |
| COD | 1.14691 | CT | 300 | 3500 | mg/l | COD | 16 mm | 3226 | 605 | OB |
| | | | 300 | 3500 | mg/l | COD | 10 mm | 4355 | | |
| COD | 1.14555 | CT | 500 | 10000 | mg/l | COD | 16 mm | 4608 | 605 | OB |
| | | | 500 | 10000 | mg/l | COD | 10 mm | 6221 | | |
| COD | 1.01797 | CT | 5000 | 90000 | mg/l | COD | 16 mm | 42553 | 605 | OB |
| | | | 5000 | 90000 | mg/l | COD | 10 mm | 57803 | | |
| COD (Hg free) | 1.09772 | CT | 10 | 150 | mg/l | COD | 16 mm | -238 | 445 | OB |
| | | | 10 | 150 | mg/l | COD | 10 mm | -321 | | |
| COD (Hg free) | 1.09773 | CT | 100 | 1500 | mg/l | COD | 16 mm | 1923 | 605 | OB |
| | | | 100 | 1500 | mg/l | COD | 10 mm | 2596 | | |
| COD for Seawater | 1.17058 | CT | 5.0 | 60.0 | mg/l | COD | 16 mm | -57.5 | 340 | OB |
| | | | 5.0 | 60.0 | mg/l | COD | 10 mm | -77.5 | | |
| COD for Seawater | 1.17059 | CT | 50 | 3000 | mg/l | COD | 16 mm | 3650 | 605 | OB |
| | | | 50 | 3000 | mg/l | COD | 10 mm | 4926 | | |
| COD | 1.18750 | CT | 0 | 40 | mg/l | COD | 16 mm | B) | 350 | OB |
| COD | 1.18751 | CT | 0 | 150 | mg/l | COD | 16 mm | B) | 420 | OB |
| COD | 1.18752 | CT | 0 | 1500 | mg/l | COD | 16 mm | B) | 620 | OB |
| COD | 1.18753 | CT | 0 | 15000 | mg/l | COD | 16 mm | B) | 620 | OB |
| Color (EN ISO 7887) | C) | | 2 | 500 | | CU | 50 mm | 380 | 410 | DW |
| | | | 0.5 | 50.0 | m ⁻¹ | | 50 mm | 20.0 | 445 | |
| | | | 0.5 | 50.0 | m ⁻¹ | | 50 mm | 20.0 | 535 | |
| | | | 0.5 | 50.0 | m ⁻¹ | | 50 mm | 20.0 | 620 | |
| Color 436 | C) | | 1 | 250 | m ⁻¹ | | 10 mm | 20.0 | 436 | DW |
| | | | 0.3 | 125.0 | m ⁻¹ | | 20 mm | 10.0 | | |
| | | | 0.1 | 50.0 | m ⁻¹ | | 50 mm | 4.00 | | |
| Color 525 | C) | | 1 | 250 | m ⁻¹ | | 10 mm | 20.0 | 525 | DW |
| | | | 0.3 | 125.0 | m ⁻¹ | | 20 mm | 10.0 | | |
| | | | 0.1 | 50.0 | m ⁻¹ | | 50 mm | 4.00 | | |
| Color 620 | C) | | 1 | 250 | m ⁻¹ | | 10 mm | 20.0 | 620 | DW |
| | | | 0.3 | 125.0 | m ⁻¹ | | 20 mm | 10.0 | | |
| | | | 0.1 | 50.0 | m ⁻¹ | | 50 mm | 4.00 | | |
| Color Measurement | see Hazen Color number | | Pt/Co or HZ | | | | | | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

B) This test can (also) be used with Hach® factory-programmed instrument calibrations. No calibration factor/table is necessary. For the corresponding Hach program number see Appendix page 18.

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaldrich.com.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|------------------|---|-----------|---------------------|-------------|------|-----------------|----------------|------------------------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Copper | 1.14767 | RT | 0.10 | 6.00 | mg/l | Cu | 10 mm | 4.44 | 605 | DW |
| | | | 0.05 | 3.00 | mg/l | Cu | 20 mm | 2.22 | | |
| | | | 0.02 | 1.20 | mg/l | Cu | 50 mm | 0.889 | | |
| | | | 0.04 | 2.36 | mg/l | Cu | 1 inch | 1.75 | | |
| Copper | 1.14553 | CT | 0.05 | 8.00 | mg/l | Cu | 16 mm | 3.60 | 605 | DW |
| | | | 0.05 | 8.00 | mg/l | Cu | 10 mm | 4.86 | | |
| Copper Bath | C) | | 10.0 | 80.0 | g/l | Cu | 10 mm | 43.1 | 820 | DW |
| | | | 5.0 | 40.0 | g/l | Cu | 20 mm | 21.6 | | |
| | | | 2.0 | 16.0 | g/l | Cu | 50 mm | 8.62 | | |
| Cyanide | 1.09701 | RT | 0.010 | 0.500 | mg/l | CN ⁻ | 10 mm | 0.221 | 605 | DW |
| | | | 0.005 | 0.2500 | mg/l | CN ⁻ | 20 mm | 0.110 | | |
| | | | 0.002 | 0.100 | mg/l | CN ⁻ | 50 mm | 0.0442 | | |
| | | | 0.004 | 0.197 | mg/l | CN ⁻ | 1 inch | 0.0871 | | |
| Cyanide | 1.14561 | CT | 0.010 | 0.500 | mg/l | CN ⁻ | 16 mm | 0.161 | 605 | DW |
| Cyanuric Acid | 1.19253 | RT | 2 | 160 | mg/l | CyA | 20 mm | B), turbidity D) | 525 | DW |
| | | | 2 | 160 | mg/l | CyA | 1 inch | | | |
| Detergents | see Surfactants | | | | | | | | | |
| Dissolved Oxygen | see Oxygen | | | | | | | | | |
| Ethanol | see Alcohol | | | | | | | | | |
| Fluoride | 1.14557 | CT | replaced by 1.00809 | | | | | | | |
| Fluoride | 1.00809 | CT | 0.10 | 1.80 | mg/l | F ⁻ | 16 mm | 1.43 | 620 | OB |
| | | | 0.10 | 1.80 | mg/l | F ⁻ | 10 mm | 1.93 | | |
| | | | 0.025 | 0.500 | mg/l | F ⁻ | 50 mm | 0.312 | | |
| Fluoride | 1.00822 | RT | 0.02 | 2.00 | mg/l | F ⁻ | 50 mm | A), B) | 605 | OB |
| | | | 0.02 | 2.00 | mg/l | F ⁻ | 1 inch | | | |
| Fluoride | 1.17236 | RT | 0.02 | 2.00 | mg/l | F ⁻ | 50 mm | A), B) | 605 | OB |
| | | | 0.02 | 2.00 | mg/l | F ⁻ | 1 inch | | | |
| Fluoride | 1.17243 | CT | 0.10 | 2.50 | mg/l | F ⁻ | 16 mm | A), B) | 605 | OB |
| | | | 0.10 | 2.50 | mg/l | F ⁻ | 50 mm | | | |
| Fluoride | 1.14598 | RT | 0.10 | 2.00 | mg/l | F ⁻ | 10 mm | 2.02 | 620 | OB |
| | | | 1.0 | 20.0 | mg/l | F ⁻ | 10 mm | 21.5 | | |
| Formaldehyde | 1.14678 | RT | 0.10 | 8.00 | mg/l | HCHO | 10 mm | 4.22 | 565 | DW |
| | | | 0.05 | 4.00 | mg/l | HCHO | 20 mm | 2.11 | | |
| | | | 0.02 | 1.50 | mg/l | HCHO | 50 mm | 0.844 | | |
| | | | 0.04 | 3.15 | mg/l | HCHO | 1 inch | 1.66 | | |
| Formaldehyde | 1.14500 | CT | 0.10 | 8.00 | mg/l | HCHO | 16 mm | 3.23 | 565 | DW |
| | | | 0.10 | 8.00 | mg/l | HCHO | 10 mm | 4.36 | | |
| Gold | 1.14821 | RT | 0.5 | 12.0 | mg/l | Au | 10 mm | 6.25 | 550 | OB |
| Hardness | see Residual Hardness or Total Hardness | | | | | | | | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

A) For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.

B) This test can (also) be used with Hach® factory-programmed instrument calibrations. No calibration factor/table is necessary. For the corresponding Hach program number see Appendix page 18.

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaaldrich.com.

D) Instrument-specific calibration needed. See Appendix page 19.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|---------------------|-----------------------|-----------|---------------------|-------------|------|-------------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Hazen Color Number | | | 0 | 500 | | CU, HZ, Pt or Pt/Co | 10 mm | 370 | 340 | DW |
| | | | 0 | 250 | | CU, HZ, Pt or Pt/Co | 20 mm | 185 | | |
| | | | 0 | 100 | | CU, HZ, Pt or Pt/Co | 50 mm | 74 | | |
| Hazen Color Number | | | 0 | 1000 | | CU, HZ, Pt or Pt/Co | 50 mm | 813 | 445 | DW |
| Hazen Color Number | | | 0 | 1000 | | CU, HZ, Pt or Pt/Co | 50 mm | 752 | 455 | DW |
| Hazen Color Number | | | 0 | 1000 | | CU, HZ, Pt or Pt/Co | 50 mm | 775 | 465 | DW |
| Hydrazine | 1.14797 | RT | replaced by 1.09711 | | | | | | | |
| Hydrazine | 1.09711 | RT | 0.02 | 2.00 | mg/l | N ₂ H ₄ | 10 mm | 0.870 | 445 | DW |
| | | | 0.01 | 1.00 | mg/l | N ₂ H ₄ | 20 mm | 0.435 | | |
| | | | 0.005 | 0.400 | mg/l | N ₂ H ₄ | 50 mm | 0.174 | | |
| | | | 0.008 | 0.787 | mg/l | N ₂ H ₄ | 1 inch | 0.342 | | |
| Hydrogen Peroxide | 1.14731 | CT | 2.0 | 20.0 | mg/l | H ₂ O ₂ | 16 mm | 38.5 | 410 | DW |
| | | | 2.0 | 20.0 | mg/l | H ₂ O ₂ | 10 mm | 52.3 | | |
| | | | 0.25 | 5.00 | mg/l | H ₂ O ₂ | 50 mm | 10.5 | | |
| Hydrogen Peroxide | 1.18789 | RT | 0.03 | 6.00 | mg/l | H ₂ O ₂ | 10 mm | 2.67 | 445 | OB |
| | | | 0.015 | 3.000 | mg/l | H ₂ O ₂ | 20 mm | 1.34 | | |
| | | | 0.01 | 2.36 | mg/l | H ₂ O ₂ | 1 inch | 1.04 | | |
| Hydrogen Sulfide | see Sulfide | | | | | | | | | |
| Hydroquinone | see Oxygen Scavengers | | | | | | | | | |
| Iodine | E) | RT | 0.20 | 10.00 | mg/l | I ₂ | 10 mm | 11.9 | 550 | DW |
| | | | 0.10 | 5.00 | mg/l | I ₂ | 20 mm | 5.95 | | |
| | | | 0.050 | 2.000 | mg/l | I ₂ | 50 mm | 2.38 | | |
| | | | 0.08 | 3.94 | mg/l | I ₂ | 1 inch | 4.69 | | |
| Iodine Color Number | | | 0.05 | 3.00 | | IFZ | 10 mm | 1.27 | 340 | DW |
| | | | 0.03 | 1.50 | | IFZ | 20 mm | 0.637 | | |
| | | | 0.010 | 0.600 | | IFZ | 50 mm | 0.255 | | |
| Iodine Color Number | | | 1.0 | 50.0 | | IFZ | 10 mm | 17.1 | 445 | DW |
| | | | 0.5 | 25.0 | | IFZ | 20 mm | 8.55 | | |
| | | | 0.2 | 10.0 | | IFZ | 50 mm | 3.42 | | |
| Iron | 1.14761 | RT | 0.05 | 5.00 | mg/l | Fe | 10 mm | 2.08 | 565 | DW |
| | | | 0.03 | 2.50 | mg/l | Fe | 20 mm | 1.04 | | |
| | | | 0.005 | 1.000 | mg/l | Fe | 50 mm | 0.416 | | |
| | | | 0.0025 | 0.5000 | mg/l | Fe | 100 mm | 0.208 | | |
| | | | 0.02 | 1.97 | mg/l | Fe | 1 inch | 0.820 | | |
| Iron | 1.00796 | RT | 0.10 | 5.00 | mg/l | Fe | 10 mm | 5.56 | 500 | DW |
| | | | 0.05 | 2.50 | mg/l | Fe | 20 mm | 2.78 | | |
| | | | 0.010 | 1.000 | mg/l | Fe | 50 mm | 1.11 | | |
| | | | 0.04 | 1.97 | mg/l | Fe | 1 inch | 2.19 | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

E) Can be determined with Chlorine test, 1.00598 (see corresponding application notes on www.sigmaaldrich.com).

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|---|-----------------------|-----------|---------------------|-------------|------|----------------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Iron | 1.14549 | CT | 0.05 | 4.00 | mg/l | Fe | 16 mm | 1.64 | 565 | DW |
| | | | 0.05 | 5.00 | mg/l | Fe | 10 mm | 2.21 | | |
| Iron | 1.14896 | CT | 1.0 | 50.0 | mg/l | Fe | 16 mm | 28.3 | 525 | DW |
| | | | 1.0 | 50.0 | mg/l | Fe | 10 mm | 38.2 | | |
| Lead | 1.09717 | RT | 0.10 | 5.00 | mg/l | Pb | 10 mm | 6.25 | 525 | OB |
| | | | 0.05 | 2.50 | mg/l | Pb | 20 mm | 3.13 | | |
| | | | 1.010 | 1.000 | mg/l | Pb | 50 mm | 1.25 | | |
| | | | 0.039 | 1.970 | mg/l | Pb | 1 inch | 2.46 | | |
| Isoascorbic Acid (Erythorbic Acid) | see Oxygen Scavengers | | | | | | | | | |
| Lead | 1.14833 | CT | 0.10 | 5.00 | mg/l | Pb | 16 mm | 4.39 | 525 | OB |
| | | | 0.10 | 5.00 | mg/l | Pb | 10 mm | 5.93 | | |
| Magnesium | 1.00815 | CT | 5.0 | 75.0 | mg/l | Mg | 16 mm | A) | 565 | DW |
| | | | 5.0 | 75.0 | mg/l | Mg | 10 mm | | | |
| Mercury | C) | | 0.025 | 1.000 | mg/l | Hg | 50 mm | 0.563 | 565 | OB |
| Manganese | 1.01739 | RT | replaced by 1.01846 | | | | | | | |
| Manganese | 1.01846 | RT | 0.05 | 2.00 | mg/l | Mn | 10 mm | 1.67 | 565 | OB |
| | | | | 1.00 | mg/l | Mn | 20 mm | 0.836 | | |
| | | | 0.005 | 0.400 | mg/l | Mn | 50 mm | 0.334 | | |
| | | | 0.020 | 0.790 | mg/l | Mn | 1 inch | 0.657 | | |
| Manganese | 1.14770 | RT | 0.50 | 10.00 | mg/l | Mn | 10 mm | 5.62 | 445 | DW |
| | | | 0.25 | 5.00 | mg/l | Mn | 20 mm | 2.81 | | |
| | | | 0.01 | 2.00 | mg/l | Mn | 50 mm | 1.12 | | |
| | | | 0.20 | 3.94 | mg/l | Mn | 1 inch | 2.21 | | |
| Manganese | 1.00816 | CT | 0.10 | 5.00 | mg/l | Mn | 16 mm | 4.13 | 445 | DW |
| | | | 0.10 | 5.00 | mg/l | Mn | 10 mm | 5.58 | | |
| Methylethylketoxime (2-Butanone-etoxime) | see Oxygen Scavengers | | | | | | | | | |
| Molybdenum | 1.00860 | CT | 0.02 | 1.00 | mg/l | Mo | 16 mm | 1.10 | 620 | OB |
| | | | 0.02 | 1.00 | mg/l | Mo | 10 mm | 1.49 | | |
| | | | 0.03 | 1.67 | mg/l | MoO ₄ ²⁻ | 16 mm | 1.84 | | |
| | | | 0.03 | 1.67 | mg/l | MoO ₄ ²⁻ | 10 mm | 2.48 | | |
| | | | 0.04 | 2.15 | mg/l | Na ₂ MoO ₄ | 16 mm | 2.37 | | |
| | | | 0.04 | 2.15 | mg/l | Na ₂ MoO ₄ | 10 mm | 3.19 | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

A) For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaldrich.com.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|----------------|-----------|-----------|------------------------------|-------------|-------|------------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Monochloramine | 1.01632 | RT | 0.25 | 10.00 | mg/l | Cl ₂ | 10 mm | 5.00 | 690 | OB |
| | | | 0.13 | 5.00 | mg/l | Cl ₂ | 20 mm | 2.50 | | |
| | | | 0.050 | 2.000 | mg/l | Cl ₂ | 50 mm | 1.00 | | |
| | | | 0.10 | 3.94 | mg/l | Cl ₂ | 1 inch | 1.97 | | |
| | | | 0.18 | 7.25 | mg/l | NH ₂ Cl | 10 mm | 3.67 | | |
| | | | 0.09 | 3.63 | mg/l | NH ₂ Cl | 20 mm | 1.83 | | |
| | | | 0.036 | 1.450 | mg/l | NH ₂ Cl | 50 mm | 0.733 | | |
| | | | 0.09 | 2.85 | mg/l | NH ₂ Cl | 1 inch | 1.44 | | |
| | | | 0.05 | 1.96 | mg/l | NH ₂ Cl-N | 10 mm | 0.988 | | |
| | | | 0.03 | 0.98 | mg/l | NH ₂ Cl-N | 20 mm | 0.494 | | |
| | | | 0.010 | 0.392 | mg/l | NH ₂ Cl-N | 50 mm | 0.198 | | |
| 0.02 | 0.77 | mg/l | NH ₂ Cl-N | 1 inch | 0.389 | | | | | |
| Nickel | 1.14785 | RT | 0.10 | 5.00 | mg/l | Ni | 10 mm | 4.76 | 445 | DW |
| | | | 0.05 | 2.50 | mg/l | Ni | 20 mm | 2.38 | | |
| | | | 0.02 | 1.00 | mg/l | Ni | 50 mm | 0.952 | | |
| | | | 0.04 | 1.97 | mg/l | Ni | 1 inch | 1.87 | | |
| Nickel | 1.14554 | CT | 0.10 | 6.00 | mg/l | Ni | 16 mm | 3.82 | 445 | OB |
| | | | 0.10 | 6.00 | mg/l | Ni | 10 mm | 5.16 | | |
| Nickel Bath | C) | | 10 | 120 | g/l | Ni | 10 mm | 60.6 | 690 | DW |
| | | | 5.0 | 60.0 | g/l | Ni | 20 mm | 30.3 | | |
| | | | 2.0 | 24.0 | g/l | Ni | 50 mm | 12.1 | | |
| Nitrate | 1.09713 | RT | 1.0 | 25.0 | mg/l | NO ₃ -N | 10 mm | 19.6 | 340 | OB |
| | | | 0.5 | 12.5 | mg/l | NO ₃ -N | 20 mm | 9.80 | | |
| | | | 0.10 | 5.00 | mg/l | NO ₃ -N | 50 mm | 3.92 | | |
| | | | 0.4 | 9.8 | mg/l | NO ₃ -N | 1 inch | 7.72 | | |
| | | | 4.4 | 110.7 | mg/l | NO ₃ ⁻ | 10 mm | 86.8 | | |
| | | | 2.2 | 55.4 | mg/l | NO ₃ ⁻ | 20 mm | 43.4 | | |
| | | | 0.4 | 22.1 | mg/l | NO ₃ ⁻ | 50 mm | 17.4 | | |
| 1.7 | 43.6 | mg/l | NO ₃ ⁻ | 1 inch | 34.2 | | | | | |
| Nitrate | 1.14773 | RT | 0.5 | 20.0 | mg/l | NO ₃ -N | 10 mm | 9.62 | 525 | OB |
| | | | 0.2 | 10.0 | mg/l | NO ₃ -N | 20 mm | 4.81 | | |
| | | | 0.2 | 7.9 | mg/l | NO ₃ -N | 1 inch | 3.79 | | |
| | | | 2.2 | 88.5 | mg/l | NO ₃ ⁻ | 10 mm | 42.6 | | |
| | | | 0.9 | 44.3 | mg/l | NO ₃ ⁻ | 20 mm | 21.3 | | |
| | | | 0.9 | 34.8 | mg/l | NO ₃ ⁻ | 1 inch | 16.8 | | |
| Nitrate | 1.14542 | CT | 0.5 | 18.0 | mg/l | NO ₃ -N | 16 mm | 7.14 | 525 | OB |
| | | | 0.5 | 18.0 | mg/l | NO ₃ -N | 10 mm | 9.64 | | |
| | | | 2.2 | 79.7 | mg/l | NO ₃ ⁻ | 16 mm | 31.6 | | |
| | | | 2.2 | 79.7 | mg/l | NO ₃ ⁻ | 10 mm | 42.7 | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaaldrich.com.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|---------------------|--------------|-----------|-----------------|-------------|------|------------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Nitrate | 1.14563 | CT | 0.5 | 25.0 | mg/l | NO ₃ -N | 16 mm | 14.7 | 340 | OB |
| | | | 0.5 | 25.0 | mg/l | NO ₃ -N | 10 mm | 19.9 | | |
| | | | 2.2 | 110.7 | mg/l | NO ₃ ⁻ | 16 mm | 65.1 | | |
| | | | 2.2 | 110.7 | mg/l | NO ₃ ⁻ | 10 mm | 87.9 | | |
| Nitrate | 1.14764 | CT | 1.0 | 50.0 | mg/l | NO ₃ -N | 16 mm | 27.9 | 340 | OB |
| | | | 1.0 | 50.0 | mg/l | NO ₃ -N | 10 mm | 37.7 | | |
| | | | 4 | 221 | mg/l | NO ₃ ⁻ | 16 mm | 124 | | |
| | | | 4 | 221 | mg/l | NO ₃ ⁻ | 10 mm | 167 | | |
| Nitrate | 1.00614 | CT | 23 | 225 | mg/l | NO ₃ -N | 16 mm | 132 | 340 | OB |
| | | | 23 | 225 | mg/l | NO ₃ -N | 10 mm | 180 | | |
| | | | 102 | 996 | mg/l | NO ₃ ⁻ | 16 mm | 586 | | |
| | | | 102 | 996 | mg/l | NO ₃ ⁻ | 10 mm | 797 | | |
| Nitrate in Seawater | 1.14556 | CT | 0.10 | 3.00 | mg/l | NO ₃ -N | 16 mm | A) | 500 | DW |
| | | | 0.10 | 2.50 | mg/l | NO ₃ -N | 10 mm | | | |
| | | | 0.4 | 13.3 | mg/l | NO ₃ ⁻ | 16 mm | | | |
| | | | 0.4 | 11.1 | mg/l | NO ₃ ⁻ | 10 mm | | | |
| Nitrate in Seawater | 1.14942 | RT | 0.2 | 17.0 | mg/l | NO ₃ ⁻ | 10 mm | A) | 500 | DW |
| | | | 0.9 | 75.3 | mg/l | NO ₃ ⁻ | 10 mm | | | |
| Nitrite | 1.14776 | RT | 0.02 | 1.00 | mg/l | NO ₂ -N | 10 mm | 0.376 | 525 | DW |
| | | | 0.010 | 0.500 | mg/l | NO ₂ -N | 20 mm | 0.188 | | |
| | | | 0.002 | 0.200 | mg/l | NO ₂ -N | 50 mm | 0.0751 | | |
| | | | 0.008 | 0.394 | mg/l | NO ₂ -N | 1 inch | 0.148 | | |
| | | | 0.07 | 3.28 | mg/l | NO ₂ ⁻ | 10 mm | 1.24 | | |
| | | | 0.03 | 1.64 | mg/l | NO ₂ ⁻ | 20 mm | 0.620 | | |
| | | | 0.007 | 0.657 | mg/l | NO ₂ ⁻ | 50 mm | 0.248 | | |
| | | | 0.026 | 1.293 | mg/l | NO ₂ ⁻ | 1 inch | 0.479 | | |
| Nitrite | 1.14547 | CT | 0.010 | 0.700 | mg/l | NO ₂ -N | 16 mm | 0.274 | 525 | DW |
| | | | 0.010 | 0.700 | mg/l | NO ₂ -N | 10 mm | 0.370 | | |
| | | | 0.03 | 2.30 | mg/l | NO ₂ -N | 16 mm | 0.900 | | |
| | | | 0.03 | 2.30 | mg/l | NO ₂ ⁻ | 10 mm | 1.22 | | |
| Nitrite | 1.00609 | CT | 1.0 | 90.0 | mg/l | NO ₂ -N | 16 mm | 82.6 | 605 | DW |
| | | | 1.0 | 90.0 | mg/l | NO ₂ -N | 10 mm | 112.4 | | |
| | | | 3.3 | 295.2 | mg/l | NO ₂ ⁻ | 16 mm | 271.3 | | |
| | | | 3.3 | 295.2 | mg/l | NO ₂ ⁻ | 10 mm | 369.2 | | |
| Nitrogen (total) | 1.00613 | CT | 0.5 | 15.0 | mg/l | N | 16 mm | 15.3 | 340 | OB |
| | | | 0.5 | 15.0 | mg/l | N | 10 mm | 20.6 | | |
| Nitrogen (total) | 1.14537 | CT | 0.5 | 15.0 | mg/l | N | 16 mm | 7.81 | 525 | OB |
| | | | 0.5 | 15.0 | mg/l | N | 10 mm | 10.5 | | |
| Nitrogen (total) | 1.14763 | CT | 10 | 150 | mg/l | N | 16 mm | 154 | 340 | OB |
| | | | 10 | 150 | mg/l | N | 10 mm | 208 | | |
| Nitrogen, Ammonium | see Ammonium | | | | | | | | | |
| Nitrogen, Nitrate | see Nitrate | | | | | | | | | |
| Nitrogen, Nitrite | see Nitrite | | | | | | | | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

A) For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|----------------------------------|----------------------------|-----------|---------------------|-------------|----------|----------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Organic Acids, Volatile | see Volatile organic acids | | | | | | | | | |
| Organic Carbon, Volatile | see TOC | | | | | | | | | |
| Oxygen, dissolved | 1.14694 | CT | 0.5 | 12.0 | mg/l | O₂ | 16 mm | 13.5 | 500 | DW |
| | | | 0.5 | 12.0 | mg/l | O₂ | 10 mm | 18.2 | | |
| Oxygen Scavengers | 1.19251 | RT | 0.020 | 0.500 | mg/l | DEHA | 20 mm | 0.408 | 565 | OB |
| | | | 0 | 450 | µg/l | DEHA | 1 inch | B) | 562 | |
| | | | 0.027 | 0.667 | mg/l | Carbohy | 20 mm | 0.544 | 565 | |
| | | | 0 | 600 | µg/l | Carbohy | 1 inch | B) | 562 | |
| | | | 0.053 | 1.315 | mg/l | Hydro | 20 mm | 1.073 | 565 | |
| | | | 0 | 1000 | µg/l | Hydro | 1 inch | B) | 562 | |
| | | | 0.078 | 1.950 | mg/l | ISA | 20 mm | 1.592 | 565 | |
| | | | 0 | 1500 | µg/l | ISA | 1 inch | B) | 562 | |
| | | | 0.087 | 2.170 | mg/l | MEKO | 20 mm | 1.771 | 565 | |
| 0 | 1000 | µg/l | MEKO | 1 inch | B) | 562 | | | | |
| Oxygen Demand, Biological | see BOD | | | | | | | | | |
| Oxygen Demand, Chemical | see COD | | | | | | | | | |
| Ozone | 1.00607 | RT | 0.05 | 4.00 | mg/l | O₃ | 10 mm | 2.25 | 550 | DW |
| | | | 0.02 | 2.00 | mg/l | O₃ | 20 mm | 1.13 | | |
| | | | 0.010 | 0.800 | mg/l | O₃ | 50 mm | 0.450 | | |
| | | | 0.020 | 1.575 | mg/l | O₃ | 1 inch | 0.887 | | |
| Ozone | 1.14732 | RT | replaced by 1.00607 | | | | | | | |
| Palladium | C) | RT | 0.05 | 1.25 | mg/l | Pd | 10 mm | 0.588 | 525 | OB |
| Peroxide | see Hydrogen Peroxide | | | | | | | | | |
| pH | 1.01744 | CT | 6.4 | 8.8 | pH units | pH | 16 mm | A) | 550 | DW |
| | | | 6.4 | 8.8 | pH units | pH | 10 mm | | | |
| Phenol | 1.00856 | RT | 0.002 | 0.100 | mg/l | Phenol | 20 mm | 0.114 | 445 | OB |
| | | | 0.10 | 5.00 | mg/l | Phenol | 10 mm | 8.33 | 500 | |
| | | | 0.002 | 0.100 | mg/l | Phenol | 1 inch | 0.0895 | 445 | |
| | | | 0.013 | 2.500 | mg/l | Phenol | 20 mm | 4.17 | 500 500 | |
| | | | 0.025 | 1.000 | mg/l | Phenol | 50 mm | 1.67 | | |
| | | | 0.07 | 1.97 | mg/l | Phenol | 1 inch | 3.28 | 495 | |
| Phenol | 1.14551 | CT | 0.10 | 2.50 | mg/l | Phenol | 16 mm | 3.39 | 500 | OB |
| | | | 0.10 | 2.50 | mg/l | Phenol | 10 mm | 4.58 | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

A) For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.

B) This test can (also) be used with Hach® factory-programmed instrument calibrations. No calibration factor/table is necessary. For the corresponding Hach program number see Appendix page 18.

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaaldrich.com.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|---------------------|-----------|-----------|-------------------------------|-------------|------|-------------------------------|----------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Phosphate (PMB) | 1.14848 | RT | 0.05 | 5.00 | mg/l | PO ₄ -P | 10 mm | 2.08 | 690 | OB |
| | | | 0.03 | 2.50 | mg/l | PO ₄ -P | 20 mm | 1.04 | | |
| | | | 0.005 | 1.000 | mg/l | PO ₄ -P | 50 mm | 0.417 | | |
| | | | 0.0025 | 0.5000 | mg/l | PO ₄ -P | 100 mm | 0.208 | | |
| | | | 0.02 | 1.97 | mg/l | PO ₄ -P | 1 inch | 0.820 | | |
| | | | 0.2 | 15.3 | mg/l | PO ₄ ³⁻ | 10 mm | 6.38 | | |
| | | | 0.09 | 7.67 | mg/l | PO ₄ ³⁻ | 20 mm | 3.19 | | |
| | | | 0.015 | 3.07 | mg/l | PO ₄ ³⁻ | 50 mm | 1.275 | | |
| | | | 0.0077 | 1.5331 | mg/l | PO ₄ ³⁻ | 100 mm | 0.638 | | |
| | | | 0.06 | 6.04 | mg/l | PO ₄ ³⁻ | 1 inch | 2.51 | | |
| | | | 0.11 | 11.46 | mg/l | P ₂ O ₅ | 10 mm | 4.76 | | |
| | | | 0.07 | 5.73 | mg/l | P ₂ O ₅ | 20 mm | 2.38 | | |
| | | | 0.02 | 2.29 | mg/l | P ₂ O ₅ | 50 mm | 0.952 | | |
| | | | 0.011 | 1.146 | mg/l | P ₂ O ₅ | 100 mm | 0.476 | | |
| 0.05 | 4.51 | mg/l | P ₂ O ₅ | 1 inch | 1.88 | | | | | |
| (o)-Phosphate (PMB) | 1.00474 | CT | 0.05 | 5.00 | mg/l | PO ₄ -P | 16 mm | 1.61 | 690 | OB |
| | | | 0.05 | 5.00 | mg/l | PO ₄ -P | 10 mm | 2.18 | | |
| | | | 0.2 | 15.3 | mg/l | PO ₄ ³⁻ | 16 mm | 4.95 | | |
| | | | 0.2 | 15.3 | mg/l | PO ₄ ³⁻ | 10 mm | 6.68 | | |
| | | | 0.11 | 11.46 | mg/l | P ₂ O ₅ | 16 mm | 3.70 | | |
| | | | 0.11 | 11.46 | mg/l | P ₂ O ₅ | 10 mm | 4.99 | | |
| Phosphate (PMB) | 1.14543 | CT | 0.05 | 5.00 | mg/l | PO ₄ -P | 16 mm | 1.61 | 690 | OB |
| | | | 0.05 | 5.00 | mg/l | PO ₄ -P | 10 mm | 2.18 | | |
| | | | 0.2 | 15.3 | mg/l | PO ₄ ³⁻ | 16 mm | 4.95 | | |
| | | | 0.2 | 15.3 | mg/l | PO ₄ ³⁻ | 10 mm | 6.68 | | |
| | | | 0.11 | 11.46 | mg/l | P ₂ O ₅ | 16 mm | 3.70 | | |
| | | | 0.11 | 11.46 | mg/l | P ₂ O ₅ | 10 mm | 4.99 | | |
| (o)-Phosphate (PMB) | 1.00475 | CT | 0.5 | 25.0 | mg/l | PO ₄ -P | 16 mm | 8.06 | 690 | OB |
| | | | 0.5 | 25.0 | mg/l | PO ₄ -P | 10 mm | 10.9 | | |
| | | | 1.5 | 76.7 | mg/l | PO ₄ ³⁻ | 16 mm | 24.7 | | |
| | | | 1.5 | 76.7 | mg/l | PO ₄ ³⁻ | 10 mm | 33.4 | | |
| | | | 1.1 | 57.3 | mg/l | P ₂ O ₅ | 16 mm | 18.5 | | |
| | | | 1.1 | 57.3 | mg/l | P ₂ O ₅ | 10 mm | 24.9 | | |
| Phosphate (PMB) | 1.14729 | CT | 0.5 | 25.0 | mg/l | PO ₄ -P | 16 mm | 8.06 | 690 | OB |
| | | | 0.5 | 25.0 | mg/l | PO ₄ -P | 10 mm | 10.9 | | |
| | | | 1.5 | 76.7 | mg/l | PO ₄ ³⁻ | 16 mm | 24.7 | | |
| | | | 1.5 | 76.7 | mg/l | PO ₄ ³⁻ | 10 mm | 33.4 | | |
| | | | 1.1 | 57.3 | mg/l | P ₂ O ₅ | 16 mm | 18.5 | | |
| | | | 1.1 | 57.3 | mg/l | P ₂ O ₅ | 10 mm | 24.9 | | |

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|--------------------------|------------------------|-----------|-----------------|-------------|------|-------------------------------|----------------|--------------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Phosphate (VM) | 1.14546 | CT | 0.5 | 25.0 | mg/l | PO ₄ -P | 16 mm | 13.4 | 410 | OB |
| | | | 0.5 | 25.0 | mg/l | PO ₄ -P | 10 mm | 18.2 | | |
| | | | 1.5 | 76.7 | mg/l | PO ₄ ³⁻ | 16 mm | 41.2 | | |
| | | | 1.5 | 76.7 | mg/l | PO ₄ ³⁻ | 10 mm | 55.6 | | |
| | | | 1.1 | 57.3 | mg/l | P ₂ O ₅ | 16 mm | 30.8 | | |
| | | | 1.1 | 57.3 | mg/l | P ₂ O ₅ | 10 mm | 41.5 | | |
| Phosphate (VM) | 1.14842 | RT | 1.0 | 30.0 | mg/l | PO ₄ -P | 10 mm | 18.0 | 410 | OB |
| | | | 0.5 | 15.0 | mg/l | PO ₄ -P | 20 mm | 9.00 | | |
| | | | 0.4 | 11.8 | mg/l | PO ₄ -P | 1 inch | 7.09 | | |
| | | | 3.1 | 92.0 | mg/l | PO ₄ ³⁻ | 10 mm | 55.2 | | |
| | | | 1.5 | 46.0 | mg/l | PO ₄ ³⁻ | 20 mm | 27.6 | | |
| | | | 1.2 | 36.2 | mg/l | PO ₄ ³⁻ | 1 inch | 21.7 | | |
| | | | 2.3 | 68.7 | mg/l | P ₂ O ₅ | 10 mm | 41.3 | | |
| | | | 1.1 | 34.4 | mg/l | P ₂ O ₅ | 20 mm | 20.6 | | |
| Phosphate (PMB) | 1.00798 | RT | 1.0 | 100.0 | mg/l | PO ₄ -P | 10 mm | 35.1 | 690 | OB |
| | | | 0.4 | 39.4 | mg/l | PO ₄ -P | 1 inch | 13.8 | | |
| | | | 3 | 307 | mg/l | PO ₄ ³⁻ | 10 mm | 108 | | |
| | | | 1.2 | 120.9 | mg/l | PO ₄ ³⁻ | 1 inch | 42.4 | | |
| | | | 2 | 229 | mg/l | P ₂ O ₅ | 10 mm | 80.4 | | |
| | | | 1 | 90 | mg/l | P ₂ O ₅ | 1 inch | 31.6 | | |
| (o)-Phosphate (PMB) | 1.00616 | CT | 3.0 | 100.0 | mg/l | PO ₄ -P | 16 mm | 39.2 | 690 | OB |
| | | | 3.0 | 100.0 | mg/l | PO ₄ -P | 10 mm | 52.9 | | |
| | | | 9 | 307 | mg/l | PO ₄ ³⁻ | 16 mm | 120 | | |
| | | | 9 | 307 | mg/l | PO ₄ ³⁻ | 10 mm | 162 | | |
| | | | 7 | 229 | mg/l | P ₂ O ₅ | 16 mm | 89.9 | | |
| | | | 7 | 229 | mg/l | P ₂ O ₅ | 10 mm | 121 | | |
| Phosphate (PMB) | 1.00673 | CT | 3.0 | 100.0 | mg/l | PO ₄ -P | 16 mm | 39.2 | 690 | OB |
| | | | 3.0 | 100.0 | mg/l | PO ₄ -P | 10 mm | 52.9 | | |
| | | | 9 | 307 | mg/l | PO ₄ ³⁻ | 16 mm | 120 | | |
| | | | 9 | 307 | mg/l | PO ₄ ³⁻ | 10 mm | 162 | | |
| | | | 7 | 229 | mg/l | P ₂ O ₅ | 16 mm | 89.9 | | |
| | | | 7 | 229 | mg/l | P ₂ O ₅ | 10 mm | 121 | | |
| Platinum | C) | RT | 0.10 | 1.25 | mg/l | Pt | 10 mm | 2.38 | 690 | OB |
| Platinum-Cobalt Standard | see Hazen Color number | | Pt/Co or HZ | | | | | | | |
| Potassium | 1.14562 | CT | 5.0 | 50.0 | mg/l | K | 16 mm | turbidity D) | 690 | DW |
| | | | 5.0 | 50.0 | mg/l | K | 10 mm | | | |
| Potassium | 1.00615 | CT | 30 | 300 | mg/l | K | 16 mm | turbidity D) | 690 | DW |
| | | | 30 | 300 | mg/l | K | 10 mm | | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaldrich.com.

D) Instrument-specific calibration needed. See Appendix page 19.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|--|-----------|-----------|-----------------|-------------|-------|-------------------|----------------------|--------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Residual Hardness | 1.14683 | CT | 0.50 | 5.00 | mg/l | Ca | 16 mm | A) | 565 | DW |
| | | | 1.25 | 12.5 | mg/l | CaCO ₃ | 16 mm | | | |
| | | | 0.70 | 7.00 | mg/l | CaO | 16 mm | | | |
| | | | 0.070 | 0.700 | | °d | 16 mm | | | |
| | | | 0.088 | 0.875 | | °e | 16 mm | | | |
| | | | 0.125 | 1.250 | | °f | 16 mm | | | |
| | | | 0.50 | 5.00 | mg/l | Ca | 10 mm | | | |
| | | | 1.25 | 12.50 | mg/l | CaCO ₃ | 10 mm | | | |
| | | | 0.70 | 7.00 | mg/l | CaO | 10 mm | | | |
| | | | 0.070 | 0.700 | | °d | 10 mm | | | |
| | | | 0.088 | 0.875 | | °e | 10 mm | | | |
| | | | 0.125 | 1.250 | | °f | 10 mm | | | |
| Silicate | 1.01813 | RT | 0.5 | 500.0 | µg/l | SiO ₂ | 50 mm | 0.599 | 820 | OB |
| | | | 0.001 | 1.000 | mg/l | SiO ₂ | 1 inch ^{*)} | 1.179 | | |
| | | | 0.2 | 233.7 | µg/l | Si | 50 mm | 0.280 | | |
| | | | 0.0004 | 0.500 | mg/l | Si | 1 inch ^{*)} | 0.551 | | |
| | | | 0.25 | 250.00 | µg/l | SiO ₂ | 100 mm | 0.299 | | |
| 0.12 | 116.85 | µg/l | Si | 100 mm | 0.140 | | | | | |
| Silicate | 1.14794 | RT | 0.21 | 10.70 | mg/l | SiO ₂ | 10 mm | 7.98 | 665 | DW |
| | | | 0.11 | 5.35 | mg/l | SiO ₂ | 20 mm | 3.99 | | |
| | | | 0.01 | 1.60 | mg/l | SiO ₂ | 50 mm | 0.611 | 820 | |
| | | | 0.08 | 4.21 | mg/l | SiO ₂ | 1 inch | 3.15 | 665 | |
| | | | 0.10 | 5.00 | mg/l | Si | 10 mm | 3.73 | 665 | |
| | | | 0.05 | 2.50 | mg/l | Si | 20 mm | 1.87 | | |
| | | | 0.005 | 0.750 | mg/l | Si | 50 mm | 0.286 | 820 | |
| 0.04 | 1.97 | mg/l | Si | 1 inch | 1.47 | 665 | | | | |
| Silicate | 1.00857 | RT | 1.1 | 107.0 | mg/l | SiO ₂ | 10 mm | 59.8 | 410 | DW |
| | | | 0.4 | 42.1 | mg/l | SiO ₂ | 1 inch | 23.5 | | |
| | | | 11 | 1070 | mg/l | SiO ₂ | 10 mm | 602 | | |
| | | | 4 | 421 | mg/l | SiO ₂ | 1 inch | 237 | | |
| | | | 0.5 | 50 | mg/l | Si | 10 mm | 27.9 | | |
| | | | 0.2 | 19.7 | mg/l | Si | 1 inch | 11.0 | | |
| | | | 5 | 500 | mg/l | Si | 10 mm | 282 | | |
| 2.0 | 196.9 | mg/l | Si | 1 inch | 111 | | | | | |
| Silver | C) | | 0.50 | 3.00 | mg/l | Ag | 10 mm | 2.22 | 550 | OB |
| | | | 0.25 | 1.50 | mg/l | Ag | 20 mm | 1.11 | | |
| | | | 0.20 | 1.18 | mg/l | Ag | 1 inch | 0.874 | | |
| Sodium in Nutrient Solutions for Fertilization | 1.00885 | CT | 10 | 300 | mg/l | Na | 16 mm | 263 | 550 | OB |
| | | | 10 | 300 | mg/l | Na | 10 mm | 355 | | |

^{*)} The programming data for the 1 inch Pour-Thru Cell and the "normal" 1-inch cell are identical.
When using the 1-inch Pour-Thru Cell please use instead of 10 ml of sample and blank 50 ml of sample and blank.

- DW** Distilled water (without reagents, poured into the type of cell used)
- OB** Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)
- A)** For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.
- C)** No ready-to-use kit. Find detailed information in the respective application note on www.sigmaldrich.com.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|-------------------------|-----------------------------|-----------|---------------------|-------------|------|-------------------------------|----------------|--------------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Sulfate | 1.01812 | RT | 2.5 | 50.0 | mg/l | SO ₄ ²⁻ | 10 mm | turbidity D) | 445 | SB |
| | | | 1.3 | 25.0 | mg/l | SO ₄ ²⁻ | 20 mm | | | |
| | | | 0.50 | 10.00 | mg/l | SO ₄ ²⁻ | 50 mm | | | |
| | | | 1.0 | 19.7 | mg/l | SO ₄ ²⁻ | 1 inch | | | |
| Sulfate | 1.02532 | CT | 1.0 | 50.0 | mg/l | SO ₄ ²⁻ | 16 mm | turbidity D) | 445 | SB |
| | | | 1.0 | 50.0 | mg/l | SO ₄ ²⁻ | 10 mm | | | |
| Sulfate | 1.14548 | CT | 5 | 250 | mg/l | SO ₄ ²⁻ | 16 mm | turbidity D) | 525 | SB |
| | | | 5 | 250 | mg/l | SO ₄ ²⁻ | 10 mm | | | |
| Sulfate | 1.02537 | RT | 5 | 300 | mg/l | SO ₄ ²⁻ | 10 mm | turbidity D) | 525 | SB |
| Sulfate | 1.00617 | CT | 50 | 500 | mg/l | SO ₄ ²⁻ | 16 mm | turbidity D) | 525 | SB |
| | | | 50 | 500 | mg/l | SO ₄ ²⁻ | 10 mm | | | |
| Sulfate | 1.14564 | CT | 100 | 1000 | mg/l | SO ₄ ²⁻ | 16 mm | turbidity D) | 820 | SB |
| | | | 100 | 1000 | mg/l | SO ₄ ²⁻ | 10 mm | | | |
| Sulfide | 1.14779 | RT | 0.10 | 1.50 | mg/l | S ²⁻ | 10 mm | A) | 665 | DW |
| | | | 0.020 | 0.500 | mg/l | S ²⁻ | 50 mm | 0.244 | | |
| Sulfite | 1.14394 | CT | 1.0 | 20.0 | mg/l | SO ₃ ²⁻ | 16 mm | 8.77 | 410 | OB |
| | | | 1.0 | 20.0 | mg/l | SO ₃ ²⁻ | 10 mm | 11.8 | | |
| | | | 0.05 | 3.00 | mg/l | SO ₃ ²⁻ | 50 mm | 1.69 | | |
| | | | 0.8 | 16.0 | mg/l | SO ₂ | 16 mm | 7.02 | | |
| | | | 0.8 | 16.0 | mg/l | SO ₂ | 10 mm | 9.44 | | |
| | | | 0.04 | 2.40 | mg/l | SO ₂ | 50 mm | 1.35 | | |
| Sulfite | 1.01746 | RT | 1.0 | 60.0 | mg/l | SO ₃ ²⁻ | 10 mm | 29.4 | 410 | OB |
| | | | 0.8 | 48.0 | mg/l | SO ₂ | 10 mm | 23.5 | | |
| Surfactants (anionic) | 1.14697 | CT | replaced by 1.02552 | | | | | | | |
| Surfactants (anionic) | 1.02552 | CT | 0.05 | 2.00 | mg/l | SDSA | 16 mm | 1.85 | 653 | OB |
| | | | 0.06 | 2.56 | mg/l | SDBS | 16 mm | 2.368 | | |
| | | | 0.05 | 2.12 | mg/l | SDS | 16 mm | 1.96 | | |
| | | | 0.08 | 3.26 | mg/l | SDOSSA | 16 mm | 3.016 | | |
| Surfactants (cationic) | 1.01764 | CT | 0.05 | 1.50 | mg/l | CTAB | 16 mm | 1.92 | 620 | OB |
| Surfactants (non-ionic) | 1.01787 | CT | 0.10 | 7.50 | mg/l | Triton X-100 | 16 mm | 6.06 | 605 | OB |
| Suspended Solids | C) | | 25 | 750 | mg/l | SuS | 20 mm | turbidity D) | 820 | DW |
| Tensides | | | | | | | | | | |
| Tin | 1.14622 | CT | replaced by 1.17265 | | | | | | | |
| Tin | 1.17265 | CT | 0.10 | 2.50 | mg/l | Sn | 16 mm | A) | 665 | DW |
| | | | 0.10 | 2.50 | mg/l | Sn | 10 mm | | | |
| TOC | 1.14878 | CT | 5.0 | 80.0 | mg/l | TOC | 16 mm | -64.5 | 605 | OB |
| TOC | 1.14879 | CT | 50 | 800 | mg/l | TOC | 16 mm | -645 | 605 | OB |
| Total Alkalinity | see Acid capacity to pH 4.3 | | | | | | | | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

SB **Sample blank** - clear sample without reagent (turbid samples must be filtered)

A) For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.

D) Instrument-specific calibration needed. See Appendix page 19.

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Factor | Wave-length [nm] | Blank |
|-------------------------------|---|-----------|--------------------------------|-------------|------|-------------------|----------------|--------------|------------------|-------|
| | | | lower limit | upper limit | | | | | | |
| Total Hardness | 1.00961 | CT | 5 | 215 | mg/l | Ca | 16 mm | A) | 565 | DW |
| | | | 13 | 538 | mg/l | CaCO ₃ | 16 mm | | | |
| | | | 7 | 301 | mg/l | CaO | 16 mm | | | |
| | | | 0.7 | 30.1 | | °d | 16 mm | | | |
| | | | 0.9 | 37.6 | | °e | 16 mm | | | |
| | | | 1.3 | 53.8 | | °f | 16 mm | | | |
| | | | 5 | 215 | mg/l | Ca | 10 mm | | | |
| | | | 13 | 538 | mg/l | CaCO ₃ | 10 mm | | | |
| | | | 7 | 301 | mg/l | CaO | 10 mm | | | |
| | | | 0.7 | 30.1 | | °d | 10 mm | | | |
| | | | 0.9 | 37.6 | | °e | 10 mm | | | |
| | | | 1.3 | 53.8 | | °f | 10 mm | | | |
| Total Nitrogen | see Nitrogen (total) | | | | | | | | | |
| Turbidity | C) | | 1 | 100 | | FAU | 50 mm | turbidity D) | 550 | DW |
| Volatile Organic Acids | 1.01763 | | replaced by 1.01749 or 1.01809 | | | | | | | |
| Volatile Organic Acids | 1.01749 | CT | 50 | 3000 | mg/l | HOAc | 16 mm | 1841 | 500 | OB |
| | | | 50 | 3000 | mg/l | HOAc | 10 mm | 2486 | | |
| | | | 73 | 4401 | mg/l | Butyric Acid | 16 mm | 2701 | | |
| | | | 73 | 4401 | mg/l | Butyric Acid | 10 mm | 3648 | | |
| Volatile Organic Acids | 1.01809 | CT | 50 | 3000 | mg/l | HOAc | 16 mm | 1841 | 500 | OB |
| | | | 50 | 3000 | mg/l | HOAc | 10 mm | 2486 | | |
| | | | 73 | 4401 | mg/l | Butyric Acid | 16 mm | 2701 | | |
| | | | 73 | 4401 | mg/l | Butyric Acid | 10 mm | 3648 | | |
| Water Hardness | see Residual Hardness or Total Hardness | | | | | | | | | |
| Zinc | 1.00861 | CT | 0.025 | 1.000 | mg/l | Zn | 16 mm | 1.36 | 500 | OB |
| | | | 0.025 | 1.000 | mg/l | Zn | 10 mm | 1.84 | | |
| Zinc | 1.14832 | RT | 0.05 | 2.50 | mg/l | Zn | 10 mm | 1.08 | 565 | DW |
| Zinc | 1.14566 | CT | 0.20 | 5.00 | mg/l | Zn | 16 mm | 4.88 | 500 | OB |
| | | | 0.20 | 5.00 | mg/l | Zn | 10 mm | 6.59 | | |

DW Distilled water (without reagents, poured into the type of cell used)

OB Own blank (distilled water and all reagents in the same quantity as used for the measuring sample preparation. Separately prepared in standard test tubes and poured into the type of cell used)

A) For programming data (non-linear calibration and absorbance-concentration-table) see Appendix page 18 and from page 20.

C) No ready-to-use kit. Find detailed information in the respective application note on www.sigmaldrich.com.

D) Instrument-specific calibration needed. See Appendix page 19.

Appendix

A) Non-linear Calibration and Absorbance-Concentration-Table

Calibration curves for the methods are non-linear.

1. Use the mentioned concentration-absorbance-tables (from page 20)
2. Besides using the concentration-absorbance-table you can enter a non-linear curve by using the coefficients of the following formula.

$$C = a + bx + cx^2 + dx^3$$

Check your instrument and enter the coefficients according the given list. The calibration is always done against distilled water.

We recommend taking a standard and checking if the calibration is done correctly because also spectrophotometers have small differences from instrument to instrument.

B) Use with Hach® factory-programmed instrument calibrations

The following tests can be used with Hach® factory-programmed instrument calibrations, so no calibration table is necessary. The corresponding Hach program number can be found in the following table:

| Parameter | Order No. | Hach program No. |
|--|-----------|------------------|
| COD | 1.18750 | 431 |
| COD | 1.18751 | 430 |
| COD | 1.18752 | 435 |
| COD | 1.18753 | 435 |
| Chlorine | 1.19254 | 80 |
| Cyanuric Acid | 1.19253 | 170 |
| Fluoride | 1.00822 | 190 |
| Fluoride | 1.17236 | 190 |
| Fluoride | 1.17243 | 190 |
| Oxygen Scavengers Carbohyrazide | 1.19251 | 180 |
| Oxygen Scavengers DEHA | 1.19251 | 181 |
| Oxygen Scavengers Hydroquinone | 1.19251 | 182 |
| Oxygen Scavengers Isoascorbic acid | 1.19251 | 183 |
| Oxygen Scavengers Methylethylketoxime | 1.19251 | 184 |

C) Alternative Methods

There are no test kits available for this parameter. The methods used here are reported in analytical literature. Applications and instructions how to perform the methods are available on the Internet homepage of Merck KGaA, Darmstadt, Germany: www.sigmaaldrich.com.

D) Turbidity Measurement

The following parameters are methods where light attenuation by turbidity is measured. As the optical system of every spectrophotometer is different, we are not proposing any factor for your spectrophotometer but suggest you prepare standards of different concentrations and determine the individual calibration factor for your respective setup.

| Parameter | Order No. |
|------------------|---------------------------------------|
| Cyanuric Acid | 1.19253 |
| Potassium | 1.14562 |
| Potassium | 1.00615 |
| Sulfate | 1.14548 |
| Sulfate | 1.00617 |
| Sulfate | 1.14564 |
| Sulfate | 1.01812 |
| Sulfate | 1.02532 |
| Sulfate | 1.02537 |
| Suspended Solids | physical measurement - application |
| Turbidity | physical measurement - application |

Acid Capacity Cell Test (Ord. No. 1.01758)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-------------------------|-----------|-----------|-----------------|-------------|--------|-------------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Acid Capacity to pH 4.3 | 1.01758 | CT | 0.40 | 8.00 | mmol/l | OH ⁻ | 16 mm | 605 | DW |
| | | | 0.40 | 8.00 | mmol/l | OH ⁻ | 10 mm | | |
| | | | 20 | 400 | mg/l | CaCO ₃ | 16 mm | | |
| | | | 20 | 400 | mg/l | CaCO ₃ | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mmol/l OH ⁻ | mg/l CaCO ₃ |
|------------------------------------|------------------------|------------------------|
| 0.462 | 0 | 0 |
| 0.488 | 0.40 | 20 |
| 0.529 | 1.00 | 50 |
| 0.593 | 1.90 | 95 |
| 0.660 | 2.80 | 140 |
| 0.728 | 3.70 | 185 |
| 0.789 | 4.50 | 225 |
| 0.856 | 5.40 | 270 |
| 0.921 | 6.30 | 315 |
| 0.984 | 7.20 | 360 |
| 1.036 | 8.00 | 400 |

| Absorbance (A) 10-mm Cell | mmol/l OH ⁻ | mg/l CaCO ₃ |
|------------------------------|------------------------|------------------------|
| 0.345 | 0 | 0 |
| 0.364 | 0.40 | 20 |
| 0.395 | 1.00 | 50 |
| 0.443 | 1.90 | 95 |
| 0.493 | 2.80 | 140 |
| 0.543 | 3.70 | 185 |
| 0.589 | 4.50 | 225 |
| 0.639 | 5.40 | 270 |
| 0.687 | 6.30 | 315 |
| 0.734 | 7.20 | 360 |
| 0.773 | 8.00 | 400 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 16-mm Round Cell | 10-mm Cell | 10-mm Cell |
|------------------------|------------------------------------|---------------------------------|------------------------------------|---------------------------------|
| | 0.40 – 8.00 mmol/l OH ⁻ | 20 – 400 mg/l CaCO ₃ | 0.40 – 8.00 mmol/l OH ⁻ | 20 – 400 mg/l CaCO ₃ |
| coefficient a: | -10.07 | -503.5 | -10.07 | -503.5 |
| coefficient b: | 30.24 | 1512 | 40.52 | 2026 |
| coefficient c: | -23.08 | -1154 | -41.44 | -2072 |
| coefficient d: | 10.36 | 517.9 | 24.93 | 1246 |

Aluminium Test (Ord. No. 1.14825)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|---------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Aluminium | 1.14825 | RT | 0.10 | 1.20 | mg/l | Al | 10 mm | 550 | DW |
| | | | 0.020 | 0.200 | mg/l | Al | 50 mm | | |
| | | | 0.039 | 0.472 | mg/l | Al | 1 inch | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 10-mm Cell | mg/l Al |
|---------------------------|---------|
| 0.189 | 0 |
| 0.248 | 0.05 |
| 0.310 | 0.10 |
| 0.374 | 0.15 |
| 0.441 | 0.20 |
| 0.585 | 0.30 |
| 0.745 | 0.40 |
| 1.110 | 0.60 |
| 1.525 | 0.80 |
| 1.936 | 1.00 |
| 2.293 | 1.20 |

| Absorbance (A) 50-mm Cell | mg/l Al |
|---------------------------|---------|
| 0.950 | 0 |
| 1.062 | 0.020 |
| 1.181 | 0.040 |
| 1.301 | 0.060 |
| 1.424 | 0.080 |
| 1.549 | 0.100 |
| 1.676 | 0.120 |
| 1.805 | 0.140 |
| 1.936 | 0.160 |
| 2.070 | 0.180 |
| 2.206 | 0.200 |

| Absorbance (A) 1-inch Cell | mg/l Al |
|----------------------------|---------|
| 0.565 | 0.040 |
| 0.680 | 0.080 |
| 0.800 | 0.120 |
| 0.958 | 0.170 |
| 1.088 | 0.210 |
| 1.259 | 0.260 |
| 1.400 | 0.300 |
| 1.547 | 0.340 |
| 1.738 | 0.390 |
| 1.901 | 0.430 |
| 2.196 | 0.500 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 10-mm Cell | 50-mm Cell | 1-inch Cell |
|------------------------|---------------------|-----------------------|-----------------------|
| | 0.10 – 1.20 mg/l Al | 0.020 – 0.240 mg/l Al | 0.040 – 0.500 mg/l Al |
| Coefficient a: | -0.1747 | -0.1747 | -0.1747 |
| Coefficient b: | 0.9866 | 0.1973 | 0.4128 |
| Coefficient c: | -0.3455 | -0.01382 | -0.06049 |
| Coefficient d: | 0.0771 | 0.000617 | 0.00565 |

Aluminium Cell Test (Ord. No. 1.00594)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|---------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Aluminium | 1.00594 | CT | 0.02 | 0.50 | mg/l | Al | 16 mm | 545 | DW |
| | | | 0.02 | 0.50 | mg/l | Al | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mg/l Al |
|------------------------------------|---------|
| 0.294 | 0 |
| 0.366 | 0.05 |
| 0.446 | 0.10 |
| 0.533 | 0.15 |
| 0.630 | 0.20 |
| 0.737 | 0.25 |
| 0.854 | 0.30 |
| 0.980 | 0.35 |
| 1.109 | 0.40 |
| 1.237 | 0.45 |
| 1.357 | 0.50 |

| Absorbance (A) 10-mm Cell | mg/l Al |
|------------------------------|---------|
| 0.219 | 0 |
| 0.273 | 0.05 |
| 0.332 | 0.10 |
| 0.398 | 0.15 |
| 0.470 | 0.20 |
| 0.550 | 0.25 |
| 0.637 | 0.30 |
| 0.731 | 0.35 |
| 0.828 | 0.40 |
| 0.923 | 0.45 |
| 1.012 | 0.50 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 10-mm Cell |
|------------------------|---------------------|---------------------|
| | 0.02 – 0.50 mg/l Al | 0.02 – 0.50 mg/l Al |
| Coefficient a: | -0.2533 | -0.2533 |
| Coefficient b: | 1.022 | 1.370 |
| Coefficient c: | -0.5990 | -1.076 |
| Coefficient d: | 0.1876 | 0.4514 |

Calcium Cell Test (Ord. No. 1.00858)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|-------------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Calcium | 1.00858 | CT | 10 | 250 | mg/l | Ca | 16 mm | 565 | DW |
| | | | 25 | 625 | mg/l | CaCO ₃ | 16 mm | | |
| | | | 14 | 350 | mg/l | CaO | 16 mm | | |
| | | | 10 | 250 | mg/l | Ca | 10 mm | | |
| | | | 25 | 625 | mg/l | CaCO ₃ | 10 mm | | |
| | | | | 350 | mg/l | CaO | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mg/l Ca | mg/l CaCO ₃ | mg/l CaO |
|------------------------------------|---------|------------------------|----------|
| 0.073 | 0 | 0 | 0 |
| 0.100 | 10 | 25 | 14 |
| 0.171 | 35 | 88 | 49 |
| 0.248 | 60 | 150 | 84 |
| 0.333 | 85 | 213 | 119 |
| 0.446 | 115 | 288 | 161 |
| 0.573 | 145 | 363 | 203 |
| 0.711 | 175 | 438 | 245 |
| 0.833 | 200 | 500 | 280 |
| 0.956 | 225 | 563 | 315 |
| 1.074 | 250 | 625 | 350 |

| Absorbance (A) 10-mm Cell | mg/l Ca | mg/l CaCO ₃ | mg/l CaO |
|------------------------------|---------|------------------------|----------|
| 0.054 | 0 | 0 | 0 |
| 0.075 | 10 | 25 | 14 |
| 0.128 | 35 | 88 | 49 |
| 0.185 | 60 | 150 | 84 |
| 0.249 | 85 | 213 | 119 |
| 0.333 | 115 | 288 | 161 |
| 0.428 | 145 | 363 | 203 |
| 0.531 | 175 | 438 | 245 |
| 0.622 | 200 | 500 | 280 |
| 0.713 | 225 | 563 | 315 |
| 0.801 | 250 | 625 | 350 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell |
|------------------------|------------------|---------------------------------|-------------------|
| | 10 – 250 mg/l Ca | 25 – 625 mg/l CaCO ₃ | 14 – 350 mg/l CaO |
| coefficient a: | -29.29 | -73.15 | -40.99 |
| coefficient b: | 416.8 | 1.041 | 583.2 |
| coefficient c: | -252.3 | -630.0 | -353.0 |
| coefficient d: | 99.16 | 247.6 | 138.7 |

| Non-linear Calibration | 10-mm Cell | 10-mm Cell | 10-mm Cell |
|------------------------|------------------|---------------------------------|-------------------|
| | 10 – 250 mg/l Ca | 25 – 625 mg/l CaCO ₃ | 14 – 350 mg/l CaO |
| coefficient a: | -29.29 | -73.15 | -40.99 |
| coefficient b: | 558.6 | 1.395 | 781.5 |
| coefficient c: | -453.0 | -1.131 | -633.8 |
| coefficient d: | 238.6 | 595.8 | 333.8 |

Chloride Test (Ord. No. 1.01807)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|-----------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Chloride | 1.01807 | RT | 0.10 | 5.00 | mg/l | Cl ⁻ | 50 mm | 500 | OB |

Absorbance-Concentration-Table, calibration against own blank

| Absorbance (A) 50-mm Cell | mg/l Cl ⁻ |
|------------------------------|----------------------|
| 0.000 | 0 |
| 0.020 | 0.10 |
| 0.062 | 0.30 |
| 0.193 | 0.90 |
| 0.329 | 1.50 |
| 0.469 | 2.10 |
| 0.607 | 2.70 |
| 0.740 | 3.30 |
| 0.867 | 3.90 |
| 0.984 | 4.50 |
| 1.075 | 5.00 |

Non-linear curve, calibration against own blank

| Non-linear Calibration | 50-mm Cell |
|------------------------|--------------------------------|
| | 0.10-5.00 mg/l Cl ⁻ |
| coefficient a: | 0.004689 |
| coefficient b: | 4.873 |
| coefficient c: | -1.357 |
| coefficient d: | 1.066 |

Chloride Cell Test (Ord. No. 1.01804)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|-----------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Chloride | 1.01804 | CT | 0.5 | 15.0 | mg/l | Cl ⁻ | 16 mm | 445 | DW |
| | | | 0.5 | 15.0 | mg/l | Cl ⁻ | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mg/l Cl ⁻ |
|------------------------------------|----------------------|
| 0.120 | 0 |
| 0.156 | 0.5 |
| 0.272 | 2.0 |
| 0.387 | 3.5 |
| 0.499 | 5.0 |
| 0.607 | 6.5 |
| 0.709 | 8.0 |
| 0.805 | 9.5 |
| 0.895 | 11.0 |
| 1.007 | 13.0 |
| 1.109 | 15.0 |

| Absorbance (A) 10-mm Cell | mg/l Cl ⁻ |
|------------------------------|----------------------|
| 0.090 | 0 |
| 0.116 | 0.5 |
| 0.203 | 2.0 |
| 0.289 | 3.5 |
| 0.373 | 5.0 |
| 0.453 | 6.5 |
| 0.529 | 8.0 |
| 0.601 | 9.5 |
| 0.668 | 11.0 |
| 0.751 | 13.0 |
| 0.828 | 15.0 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 10-mm Cell |
|------------------------|-------------------------------|-------------------------------|
| | 0.5-15.0 mg/l Cl ⁻ | 0.5-15.0 mg/l Cl ⁻ |
| coefficient a: | -1.537 | -1.537 |
| coefficient b: | 13.35 | 17.89 |
| coefficient c: | -2.082 | -3.738 |
| coefficient d: | 3.146 | 7.570 |

Fluoride Test (Ord. No. 1.00822)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|----------------|----------------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Fluoride | 1.00822 | RT | 0.02 | 2.00 | mg/l | F ⁻ | 50 mm | 605 | OB |
| | | | 0.02 | 2.00 | mg/l | F ⁻ | 1 inch ^{*)} | | |

Absorbance-Concentration-Table, calibration against own blank

| Absorbance (A) 50-mm Cell | mg/l F ⁻ |
|------------------------------|---------------------|
| 0.001 | 0 |
| -0.010 | 0.02 |
| -0.146 | 0.25 |
| -0.296 | 0.50 |
| -0.447 | 0.75 |
| -0.594 | 1.00 |
| -0.708 | 1.20 |
| -0.816 | 1.40 |
| -0.920 | 1.60 |
| -1.018 | 1.80 |
| -1.111 | 2.00 |

Non-linear curve, calibration against own blank

| Non-linear Calibration | 50-mm Cell |
|------------------------|---------------------------------|
| | 0.02 – 2.00 mg/l F ⁻ |
| coefficient a: | 0.002 |
| coefficient b: | -1.729 |
| coefficient c: | -0.253 |
| coefficient d: | -0.284 |

*) The Fluoride Test 1.00822 can be used with Hach factory-programmed instrument calibrations (Hach program no. 190). No calibration table is necessary.

Handling is identical to the corresponding Hach test. Details can be found in the package insert, which is included in the box.

Fluoride Test (Ord. No. 1.17236)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|----------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Fluoride | 1.17236 | RT | 0.02 | 2.00 | mg/l | F ⁻ | 50 mm | 605 | OB |
| | | | 0.02 | 2.00 | mg/l | F ⁻ | 1 inch | | |

Absorbance-Concentration-Table, calibration against own blank

| Absorbance (A) 50-mm Cell | mg/l F ⁻ |
|------------------------------|---------------------|
| 0.000 | 0.00 |
| -0.018 | 0.02 |
| -0.155 | 0.25 |
| -0.307 | 0.50 |
| -0.458 | 0.75 |
| -0.605 | 1.00 |
| -0.730 | 1.20 |
| -0.842 | 1.40 |
| -0.950 | 1.60 |
| -1.054 | 1.80 |
| -1.141 | 2.00 |

Non-linear curve, calibration against own blank

| Non-linear Calibration | 50-mm Cell |
|------------------------|---------------------------------|
| | 0.02 – 2.00 mg/l F ⁻ |
| coefficient a: | 0.002385 |
| coefficient b: | -1.729 |
| coefficient c: | -0.2529 |
| coefficient d: | -0.2843 |

*) The Fluoride Test 1.17236 can be used with Hach factory-programmed instrument calibrations (Hach program no. 190). No calibration table is necessary.

Handling is identical to the corresponding Hach test. Details can be found in the package insert, which is included in the box.

Fluoride Cell Test (Ord. No. 1.17243)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|----------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Fluoride | 1.17243 | CT | 0.10 | 2.50 | mg/l | F ⁻ | 16 mm | 605 | OB |
| | | | 0.10 | 2.50 | mg/l | F ⁻ | 50 mm | | |

Absorbance-Concentration-Table, calibration against own blank

| Absorbance (A) 16-mm Round Cell | mg/l F ⁻ |
|---------------------------------------|---------------------|
| 0.000 | 0.00 |
| -0.030 | 0.10 |
| -0.072 | 0.30 |
| -0.117 | 0.50 |
| -0.168 | 0.75 |
| -0.225 | 1.00 |
| -0.288 | 1.30 |
| -0.345 | 1.60 |
| -0.410 | 2.00 |
| -0.441 | 2.25 |
| -0.464 | 2.50 |

| Absorbance (A) 50-mm Cell | mg/l F ⁻ |
|---------------------------------|---------------------|
| 0.000 | 0.00 |
| -0.111 | 0.10 |
| -0.268 | 0.30 |
| -0.434 | 0.50 |
| -0.623 | 0.75 |
| -0.832 | 1.00 |
| -1.065 | 1.30 |
| -1.276 | 1.60 |
| -1.519 | 2.00 |
| -1.635 | 2.25 |
| -1.719 | 2.50 |

Non-linear curve, calibration against own blank

| Non-linear Calibration | 16-mm Round Cell | 50-mm Cell |
|---------------------------|------------------------------------|------------------------------------|
| | 0.10 – 2.50 mg/l F ⁻ | 0.10 – 2.50 mg/l F ⁻ |
| coefficient a: | -0.02737 | -0.02737 |
| coefficient b: | -5.140 | -1.388 |
| coefficient c: | -6.015 | -0.4385 |
| coefficient d: | -14.00 | -0.2755 |

*) The Fluoride Cell Test 1.17243 can be used with Hach factory-programmed instrument calibrations (Hach program no. 190). No calibration table is necessary.

Handling is identical to the corresponding Hach test. Details can be found in the package insert, which is included in the box.

Magnesium Cell Test (Ord. No. 1.00815)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|---------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Magnesium | 1.00815 | CT | 5.0 | 75.0 | mg/l | Mg | 16 mm | 565 | DW |
| | | | 5.0 | 75.0 | mg/l | Mg | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mg/l Mg |
|---------------------------------------|---------|
| 0.069 | 0 |
| 0.087 | 5.0 |
| 0.119 | 13.0 |
| 0.151 | 20.0 |
| 0.190 | 28.0 |
| 0.235 | 36.0 |
| 0.284 | 44.0 |
| 0.337 | 52.0 |
| 0.384 | 59.0 |
| 0.436 | 67.0 |
| 0.484 | 75.0 |

| Absorbance (A) 10-mm Cell | mg/l Mg |
|---------------------------------|---------|
| 0.051 | 0 |
| 0.065 | 5.0 |
| 0.089 | 13.0 |
| 0.112 | 20.0 |
| 0.142 | 28.0 |
| 0.175 | 36.0 |
| 0.212 | 44.0 |
| 0.251 | 52.0 |
| 0.287 | 59.0 |
| 0.325 | 67.0 |
| 0.361 | 75.0 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 10-mm Cell |
|------------------------|--------------------|--------------------|
| | 5.0 – 75.0 mg/l Mg | 5.0 – 75.0 mg/l Mg |
| coefficient a: | -21.51 | -21.5 |
| coefficient b: | 350.6 | 469.7 |
| coefficient c: | -577.9 | -1038 |
| coefficient d: | 548.8 | 1320 |

Nitrate Cell Test in seawater (Ord. No. 1.14556)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|---------------------|-----------|-----------|-----------------|-------------|------|-----------------------------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Nitrate in Seawater | 1.14556 | CT | 0.10 | 3.00 | mg/l | NO₃-N | 16 mm | 500 | DW |
| | | | 0.10 | 2.50 | mg/l | NO₃-N | 10 mm | | |
| | | | 0.4 | 13.3 | mg/l | NO₃⁻ | 16 mm | | |
| | | | 0.4 | 11.1 | mg/l | NO₃⁻ | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mg/l NO ₃ -N | mg/l NO ₃ ⁻ |
|------------------------------------|-------------------------|-----------------------------------|
| 0.000 | 0 | 0 |
| 0.088 | 0.10 | 0.4 |
| 0.118 | 0.15 | 0.7 |
| 0.279 | 0.40 | 1.8 |
| 0.492 | 0.70 | 3.1 |
| 0.726 | 1.00 | 4.4 |
| 0.986 | 1.30 | 5.8 |
| 1.275 | 1.60 | 7.1 |
| 1.583 | 1.90 | 8.4 |
| 1.904 | 2.20 | 9.7 |
| 2.216 | 2.50 | 11.1 |
| 2.510 | 2.80 | 12.4 |
| 2.690 | 3.00 | 13.3 |

| Absorbance (A) 10-mm Cell | mg/l NO ₃ -N | mg/l NO ₃ ⁻ |
|------------------------------|-------------------------|-----------------------------------|
| 0.000 | 0 | 0 |
| 0.066 | 0.10 | 0.4 |
| 0.088 | 0.15 | 0.7 |
| 0.208 | 0.40 | 1.8 |
| 0.367 | 0.70 | 3.1 |
| 0.542 | 1.00 | 4.4 |
| 0.736 | 1.30 | 5.8 |
| 0.951 | 1.60 | 7.1 |
| 1.181 | 1.90 | 8.4 |
| 1.421 | 2.20 | 9.7 |
| 1.654 | 2.50 | 11.1 |
| 1.873 | 2.80 | 12.4 |
| 2.007 | 3.00 | 13.3 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 16-mm Round Cell | 10-mm Cell | 10-mm Cell |
|------------------------|-------------------------------------|--|-------------------------------------|--|
| | 0.10 – 3.00 mg/l NO ₃ -N | 0.4 – 13.3 mg/l NO ₃ ⁻ | 0.10 – 3.00 mg/l NO ₃ -N | 0.4 – 13.3 mg/l NO ₃ ⁻ |
| coefficient a: | -0.0476 | -0.2107 | -0.04760 | -0.2107 |
| coefficient b: | 1.718 | 7.604 | 2.302 | 10.19 |
| coefficient c: | -0.4382 | -1.940 | -0.7868 | -3.483 |
| coefficient d: | 0.08210 | 0.3634 | 0.1975 | 0.8745 |

Nitrate Test in seawater (Ord. No. 1.14942)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|---------------------|-----------|-----------|-----------------|-------------|------|------------------------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Nitrate in Seawater | 1.14942 | RT | 0.2 | 17.0 | mg/l | NO ₃ ⁻ | 10 mm | 500 | DW |
| | | | 0.9 | 75.3 | mg/l | NO ₃ ⁻ | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 10-mm Cell | mg/l NO ₃ -N | mg/l NO ₃ ⁻ |
|---------------------------|-------------------------|-----------------------------------|
| 0.000 | 0.0 | 0.0 |
| 0.070 | 1.0 | 4.4 |
| 0.195 | 2.0 | 8.9 |
| 0.320 | 3.0 | 13.3 |
| 0.575 | 5.0 | 22.1 |
| 0.845 | 7.0 | 31.0 |
| 1.120 | 9.0 | 39.8 |
| 1.407 | 11.0 | 48.7 |
| 1.706 | 13.0 | 57.6 |
| 2.012 | 15.0 | 66.4 |
| 2.335 | 17.0 | 75.3 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 10-mm Cell | 10-mm Cell |
|------------------------|------------------------------------|--|
| | 0.2 – 17.0 mg/l NO ₃ -N | 0.9 – 75.3 mg/l NO ₃ ⁻ |
| coefficient a: | 0.4213 | 1.865 |
| coefficient b: | 8.251 | 36.53 |
| coefficient c: | -0.5679 | -2.514 |
| coefficient d: | 0.0335 | 0.1481 |

pH Cell Test (Ord. No. 1.01744)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|----------|---------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| pH | 1.01744 | CT | 6.4 | 8.8 | pH units | pH | 16 mm | 550 | DW |
| | | | 6.4 | 8.8 | pH units | pH | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | pH |
|------------------------------------|-----|
| 0.000 | 6.2 |
| 0.068 | 6.4 |
| 0.192 | 6.7 |
| 0.345 | 7.0 |
| 0.470 | 7.2 |
| 0.615 | 7.4 |
| 0.800 | 7.6 |
| 1.020 | 7.8 |
| 1.260 | 8.0 |
| 1.640 | 8.4 |
| 1.890 | 8.8 |

| Absorbance (A) 10-mm Cell | pH |
|------------------------------|-----|
| 0.000 | 6.2 |
| 0.051 | 6.4 |
| 0.143 | 6.7 |
| 0.257 | 7.0 |
| 0.351 | 7.2 |
| 0.459 | 7.4 |
| 0.597 | 7.6 |
| 0.761 | 7.8 |
| 0.940 | 8.0 |
| 1.224 | 8.4 |
| 1.410 | 8.8 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 10-mm Cell |
|------------------------|------------------|--------------|
| | pH 6.4 – 8.8 | pH 6.4 – 8.8 |
| coefficient a: | 6.217 | 6.217 |
| coefficient b: | 2.836 | 3.800 |
| coefficient c: | -1.824 | -3.275 |
| coefficient d: | 0.5531 | 1.331 |

Residual Hardness Cell Test (Ord. No. 1.14683)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-------------------|-----------|-----------|-----------------|-------------|------|-------------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Residual Hardness | 1.14683 | CT | 0.50 | 5.00 | mg/l | Ca | 16 mm | 565 | DW |
| | | | 1.25 | 12.5 | mg/l | CaCO ₃ | 16 mm | | |
| | | | 0.70 | 7.00 | mg/l | CaO | 16 mm | | |
| | | | 0.070 | 0.700 | | °d | 16 mm | | |
| | | | 0.088 | 0.875 | | °e | 16 mm | | |
| | | | 0.125 | 1.250 | | °f | 16 mm | | |
| | | | 0.50 | 5.00 | mg/l | Ca | 10 mm | | |
| | | | 1.25 | 12.50 | mg/l | CaCO ₃ | 10 mm | | |
| | | | 0.70 | 7.00 | mg/l | CaO | 10 mm | | |
| | | | 0.070 | 0.700 | | °d | 10 mm | | |
| 0.088 | 0.875 | | °e | 10 mm | | | | | |
| 0.125 | 1.250 | | °f | 10 mm | | | | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mg/l Ca | mg/l CaCO ₃ | mg/l CaO | °d | °e | °f |
|---------------------------------------|---------|------------------------|----------|-------|-------|------|
| 0.100 | 0 | 0.0 | 0 | 0 | 0 | 0.00 |
| 0.219 | 0.50 | 1.3 | 0.70 | 0.070 | 0.088 | 0.13 |
| 0.343 | 1.00 | 2.5 | 1.40 | 0.140 | 0.175 | 0.25 |
| 0.473 | 1.50 | 3.8 | 2.10 | 0.210 | 0.263 | 0.38 |
| 0.609 | 2.00 | 5.0 | 2.80 | 0.280 | 0.350 | 0.50 |
| 0.751 | 2.50 | 6.3 | 3.50 | 0.350 | 0.438 | 0.63 |
| 0.898 | 3.00 | 7.5 | 4.20 | 0.420 | 0.525 | 0.75 |
| 1.201 | 4.00 | 10.0 | 5.60 | 0.560 | 0.700 | 1.00 |
| 1.503 | 5.00 | 12.5 | 7.00 | 0.700 | 0.875 | 1.25 |

| Absorbance (A) 10-mm Cell | mg/l Ca | mg/l CaCO ₃ | mg/l CaO | °d | °e | °f |
|---------------------------------|---------|------------------------|----------|-------|-------|------|
| 0.075 | 0 | 0.0 | 0 | 0 | 0 | 0.00 |
| 0.164 | 0.50 | 1.3 | 0.70 | 0.070 | 0.088 | 0.13 |
| 0.255 | 1.00 | 2.5 | 1.40 | 0.140 | 0.175 | 0.25 |
| 0.353 | 1.50 | 3.8 | 2.10 | 0.210 | 0.263 | 0.38 |
| 0.454 | 2.00 | 5.0 | 2.80 | 0.280 | 0.350 | 0.50 |
| 0.560 | 2.50 | 6.3 | 3.50 | 0.350 | 0.438 | 0.63 |
| 0.670 | 3.00 | 7.5 | 4.20 | 0.420 | 0.525 | 0.75 |
| 0.896 | 4.00 | 10.0 | 5.60 | 0.560 | 0.700 | 1.00 |
| 1.121 | 5.00 | 12.5 | 7.00 | 0.700 | 0.875 | 1.25 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell |
|------------------------|---------------------|-------------------------------------|----------------------|------------------|------------------|------------------|
| | 0.50 – 5.00 mg/l Ca | 1.25 – 12.50 mg/l CaCO ₃ | 0.70 – 7.00 mg/l CaO | 0.070 – 0.700 °d | 0.088 – 0.875 °e | 0.125 – 1.250 °f |
| coefficient a: | -0.4615 | -1.152 | -0.6457 | -0.06458 | -0.08065 | -0.1153 |
| coefficient b: | 4.623 | 11.54 | 6.467 | 0.6468 | 0.8079 | 1.154 |
| coefficient c: | -1.141 | -2.849 | -1.596 | -0.1596 | -0.1994 | -0.2849 |
| coefficient d: | 0.3236 | 0.8080 | 0.4527 | 0.04528 | 0.05655 | 0.08081 |

| Non-linear Calibration | 10-mm Cell | 10-mm Cell | 10-mm Cell | 10-mm Cell | 10-mm Cell | 10-mm Cell |
|------------------------|---------------------|-------------------------------------|----------------------|------------------|------------------|------------------|
| | 0.50 – 5.00 mg/l Ca | 1.25 – 12.50 mg/l CaCO ₃ | 0.70 – 7.00 mg/l CaO | 0.070 – 0.700 °d | 0.088 – 0.875 °e | 0.125 – 1.250 °f |
| coefficient a: | -0.4615 | -1.152 | -0.6457 | -0.06458 | -0.08065 | -0.1153 |
| coefficient b: | 6.195 | 15.47 | 8.666 | 0.8668 | 1.083 | 1.547 |
| coefficient c: | -2.048 | -5.115 | -2.866 | -0.2866 | -0.3580 | -0.5116 |
| coefficient d: | 0.7786 | 1.944 | 1.089 | 0.1089 | 0.1361 | 0.1944 |

Sulfide Test (Ord. No. 1.14779)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|-----------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Sulfide | 1.14779 | RT | 0.10 | 1.50 | mg/l | S ²⁻ | 10 mm | 665 | DW |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 10-mm Cell | mg/l S ²⁻ |
|------------------------------|----------------------|
| 0.000 | 0 |
| 0.078 | 0.10 |
| 0.204 | 0.25 |
| 0.328 | 0.40 |
| 0.450 | 0.55 |
| 0.569 | 0.70 |
| 0.720 | 0.90 |
| 0.827 | 1.05 |
| 1.028 | 1.35 |
| 1.121 | 1.50 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 10-mm Cell |
|------------------------|----------------------------------|
| | 0.10 – 1.50 mg/l S ²⁻ |
| coefficient a: | 0.007200 |
| coefficient b: | 1.193 |
| coefficient c: | -0.03710 |
| coefficient d: | 0.1430 |

Tin Cell Test (Ord. No. 1.17265)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|-----------|-----------|-----------|-----------------|-------------|------|---------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Tin | 1.17265 | CT | 0.10 | 2.50 | mg/l | Sn | 16 mm | 665 | DW |
| | | | 0.10 | 2.50 | mg/l | Sn | 10 mm | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mg/l Sn |
|---------------------------------------|---------|
| 0.000 | 0.00 |
| 0.067 | 0.10 |
| 0.294 | 0.40 |
| 0.529 | 0.70 |
| 0.765 | 1.00 |
| 0.937 | 1.20 |
| 1.180 | 1.50 |
| 1.361 | 1.70 |
| 1.625 | 2.00 |
| 1.860 | 2.30 |
| 1.989 | 2.50 |

| Absorbance (A) 10-mm Cell | mg/l Sn |
|---------------------------------|---------|
| 0.000 | 0.00 |
| 0.049 | 0.10 |
| 0.218 | 0.40 |
| 0.391 | 0.70 |
| 0.566 | 1.00 |
| 0.694 | 1.20 |
| 0.874 | 1.50 |
| 1.008 | 1.70 |
| 1.204 | 2.00 |
| 1.378 | 2.30 |
| 1.473 | 2.50 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 10-mm Cell |
|------------------------|---------------------|---------------------|
| | 0.10 – 2.50 mg/l Sn | 0.10 – 2.50 mg/l Sn |
| coefficient a: | -0.001857 | -0.001857 |
| coefficient b: | 1.466 | 1.979 |
| coefficient c: | -0.2730 | -0.4976 |
| coefficient d: | 0.08289 | 0.2039 |

Total Hardness Cell Test (Ord. No. 1.00961)

| Parameter | Order No. | Test Type | Measuring range | | Unit | Chemical Form | Cell size used | Wave-length [nm] | Blank |
|----------------|-----------|-----------|-----------------|-------------|------|-------------------|----------------|------------------|-------|
| | | | lower limit | upper limit | | | | | |
| Total Hardness | 1.00961 | CT | 5 | 215 | mg/l | Ca | 16 mm | 565 | DW |
| | | | 13 | 538 | mg/l | CaCO ₃ | 16 mm | | |
| | | | 7 | 301 | mg/l | CaO | 16 mm | | |
| | | | 0.7 | 30.1 | | °d | 16 mm | | |
| | | | 0.9 | 37.6 | | °e | 16 mm | | |
| | | | 1.3 | 53.8 | | °f | 16 mm | | |
| | | | 5 | 215 | mg/l | Ca | 10 mm | | |
| | | | 13 | 538 | mg/l | CaCO ₃ | 10 mm | | |
| | | | 7 | 301 | mg/l | CaO | 10 mm | | |
| | | | 0.7 | 30.1 | | °d | 10 mm | | |
| 0.9 | 37.6 | | °e | 10 mm | | | | | |
| 1.3 | 53.8 | | °f | 10 mm | | | | | |

Absorbance-Concentration-Table, calibration against distilled water

| Absorbance (A) 16-mm Round Cell | mg/l Ca | mg/l CaCO ₃ | mg/l CaO | °d | °e | °f |
|---------------------------------|---------|------------------------|----------|------|------|------|
| 0.068 | 0 | 0 | 0 | 0 | 0 | 0 |
| 0.085 | 5 | 13 | 7 | 0.7 | 0.9 | 1.3 |
| 0.152 | 25 | 63 | 35 | 3.5 | 4.4 | 6.3 |
| 0.245 | 50 | 125 | 70 | 7.0 | 8.8 | 12.5 |
| 0.348 | 75 | 188 | 105 | 10.5 | 13.1 | 18.8 |
| 0.464 | 100 | 250 | 140 | 14.0 | 17.5 | 25.0 |
| 0.567 | 120 | 300 | 168 | 16.8 | 21.0 | 30.0 |
| 0.708 | 145 | 363 | 203 | 20.3 | 25.4 | 36.3 |
| 0.862 | 170 | 425 | 238 | 23.8 | 29.8 | 42.5 |
| 1.021 | 195 | 488 | 273 | 27.3 | 34.1 | 48.8 |
| 1.144 | 215 | 538 | 301 | 30.1 | 37.6 | 53.8 |

| Absorbance (A) 10-mm Cell | mg/l Ca | mg/l CaCO ₃ | mg/l CaO | °d | °e | °f |
|---------------------------|---------|------------------------|----------|------|------|------|
| 0.051 | 0 | 0 | 0 | 0 | 0 | 0 |
| 0.063 | 5 | 13 | 7 | 0.7 | 0.9 | 1.3 |
| 0.113 | 25 | 63 | 35 | 3.5 | 4.4 | 6.3 |
| 0.183 | 50 | 125 | 70 | 7.0 | 8.8 | 12.5 |
| 0.260 | 75 | 188 | 105 | 10.5 | 13.1 | 18.8 |
| 0.346 | 100 | 250 | 140 | 14.0 | 17.5 | 25.0 |
| 0.423 | 120 | 300 | 168 | 16.8 | 21.0 | 30.0 |
| 0.528 | 145 | 363 | 203 | 20.3 | 25.4 | 36.3 |
| 0.643 | 170 | 425 | 238 | 23.8 | 29.8 | 42.5 |
| 0.762 | 195 | 488 | 273 | 27.3 | 34.1 | 48.8 |
| 0.854 | 215 | 538 | 301 | 30.1 | 37.6 | 53.8 |

Non-linear curve, calibration against distilled water

| Non-linear Calibration | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell | 16-mm Round Cell |
|------------------------|------------------|---------------------------------|------------------|------------------|------------------|------------------|
| | 5 – 215 mg/l Ca | 13 – 538 mg/l CaCO ₃ | 7 – 301 mg/l CaO | 0.7 – 30.1 °d | 0.9 – 37.6 °e | 1.3 – 53.8 °f |
| coefficient a: | -22.35 | -55.82 | -31.27 | -3.128 | -3.907 | -5.582 |
| coefficient b: | 340.0 | 848.9 | 475.6 | 47.57 | 59.41 | 84.90 |
| coefficient c: | -195.3 | -487.6 | -273.2 | -27.32 | -34.12 | -48.76 |
| coefficient d: | 69.67 | 174.0 | 97.47 | 9.748 | 12.18 | 17.40 |

| Non-linear Calibration | 10-mm Cell | 10-mm Cell | 10-mm Cell | 10-mm Cell | 10-mm Cell | 10-mm Cell |
|------------------------|-----------------|---------------------------------|------------------|---------------|---------------|---------------|
| | 5 – 215 mg/l Ca | 13 – 538 mg/l CaCO ₃ | 7 – 301 mg/l CaO | 0.7 – 30.1 °d | 0.9 – 37.6 °e | 1.3 – 53.8 °f |
| coefficient a: | -22.35 | -55.82 | -31.27 | -3.128 | -3.907 | -5.582 |
| coefficient b: | 455.6 | 1.138 | 637.3 | 63.74 | 79.61 | 113.8 |
| coefficient c: | -350.6 | -875.5 | -490.5 | -49.06 | -61.27 | -87.55 |
| coefficient d: | 167.6 | 418.6 | 234.5 | 23.46 | 29.30 | 41.86 |

Merck KGaA, 64271 Darmstadt, Germany, Tel. +49(0)6151 72-2440
www.sigmaldrich.com
EMD Millipore Corporation, 400 Summit Drive, Burlington MA 01803, USA,
Tel. +1-978-715-4321
Sigma-Aldrich Canada Co. or Millipore (Canada) Ltd. 2149 Winston Park,
Dr. Oakville, Ontario, L6H 6J8, Phone: +1 800-565-1400

MilliporeSigma is the U.S. and Canada Life Science business of Merck KGaA, Darmstadt, Germany.

© 2023 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. MilliporeSigma, the vibrant M, Supelco, Sigma-Aldrich, and Spectroquant are trademarks of Merck KGaA, Darmstadt, Germany. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly available resources.

